



Metabolism Assay Kits

Convenient Assay Kits to Analyze Metabolites and Enzymes

Features and Benefits

- Convenient, simple and highly dependable assays for monitoring metabolic pathways
- Assay kits utilize spectrophotometric, fluorometric and/or luminescence detection methods
- Most assay kits are suitable for high-throughput assays

Assay Kits for Quantification of Metabolites and Enzymes

We offer a wide range of kits for analyzing both critical metabolites and the activity of key metabolic enzymes.

Name	Description	Cat No.
Glycolysis Metabolism Assay Kits		
Acetate Colorimetric Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK474
Acetyl-Coenzyme A Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK566
ADP Assay Kit	sufficient for 100 fluorometric tests	» MAK518
ATP Assay Kit	sufficient for 100 assays (bioluminescent)	» MAK473
Cytotoxicity Assay Kit	sufficient for 100 assays (bioluminescent)	» MAK536
Glucose-6-Phosphate Dehydrogenase Assay Kit	sufficient for 100 colorimetric tests	» MAK451
Glucose Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK476
Glucose (GO) Assay Kit	sufficient for 20 tests	» GAGO20
Glucose (HK) Assay Kit	sufficient for 20 tests	» GAHK020
Glucose Uptake Assay Kit	sufficient for 100 colorimetric tests	» MAK542
Glucose Uptake Assay Kit	sufficient for 100 fluorometric tests	» MAK489
Glucose-6-Phosphate Assay Kit	sufficient for 100 colorimetric tests	» MAK503

Name	Description	Cat No.
Glucose-6-Phosphate (G6P) Assay Kit	sufficient for 200 fluorometric tests	» MAK548
Glycerol 3-Phosphate (G3P) Assay Kit	sufficient for 200 colorimetric tests	» MAK546
Glycogen Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK465
High Sensitivity Glucose Quantitation Kit	sufficient for 500 fluorometric tests	» MAK543
L-Lactate Assay Kit	sufficient for 100 colorimetric tests	» MAK329
L-Lactate Assay Kit	sufficient for 100 fluorometric tests	» MAK443
Lactate Dehydrogenase Assay Kit	sufficient for 100 colorimetric tests	» MAK464
LDH Cytotoxicity Assay Kit	sufficient for 100 colorimetric tests	» MAK529
Lactose Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK487
Malate Assay Kit	sufficient for 100 colorimetric tests	» MAK511
Oxaloacetate Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK515

Name	Description	Cat No.
Pyruvate Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK332
Pyruvate Kinase Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK485
Total Carbohydrate Assay Kit	sufficient for 100 colorimetric tests	» MAK559
TCA Metabolism Assay Kits		
Acetyl-Coenzyme A Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK566
Aconitase Assay Kit	sufficient for 100 colorimetric tests	» MAK337
Citrate Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK333
Citrate Synthase Assay Kit	sufficient for 480 colorimetric reactions	» CS0720
Coenzyme A Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK504
Fumarate Assay Kit	sufficient for 100 colorimetric tests	» MAK492
Isocitrate Dehydrogenase (IDH) Assay Kit	sufficient for 100 colorimetric tests	» MAK493
Isocitrate Assay Kit	sufficient for 100 colorimetric tests	» MAK524
Malate Assay Kit	sufficient for 100 colorimetric tests	» MAK511
Malate Dehydrogenase Assay Kit	sufficient for 100 colorimetric tests	» MAK512
Oxaloacetate Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK515
α -Ketoglutarate Quantitation Kit	sufficient for 200 colorimetric tests	» MAK541
Oxidative Stress Metabolism Assay Kits		
Ascorbic Acid Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK505
Carbonyl Assay Kit	sufficient for 100 colorimetric tests	» MAK486
Fluorometric Intracellular ROS KIT	sufficient for 200 fluorometric tests (deep red)	» MAK142
Fluorometric Intracellular ROS KIT	sufficient for 200 fluorometric tests (green)	» MAK143
Fluorometric Intracellular ROS KIT	sufficient for 200 fluorometric tests (orange)	» MAK144
Fluorometric Intracellular ROS KIT	sufficient for 200 fluorometric tests (red)	» MAK145
Hemoglobin Assay Kit	sufficient for 250 colorimetric tests	» MAK115
Hypochlorite (Hypochlorous Acid) Assay Kit	sufficient for 200 colorimetric tests	» MAK556
Lipid Peroxidation (MDA) Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK568
Malate Dehydrogenase Assay Kit	sufficient for 100 colorimetric tests	» MAK512

Name	Description	Cat No.
Melanin Assay Kit	sufficient for 100 fluorometric tests	» MAK557
Monoamine Oxidase Activity Assay Kit	sufficient for 100 fluorometric tests	» MAK136
MPO Colorimetric Activity Assay Kit	sufficient for 100 colorimetric tests	» MAK563
MPO Fluorometric Activity Assay Kit	sufficient for 100 fluorometric tests	» MAK516
Nitrite/Nitrate Assay Kit, colorimetric	sufficient for 100 colorimetric tests	» 23479
Peroxidase Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK506
Superoxide Dismutase (SOD) Assay Kit	sufficient for 100 colorimetric tests	» MAK528
Uric Acid Assay Kit	sufficient for 250 colorimetric tests	» MAK483
Xanthine Oxidase Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK497
Cholesterol Metabolism Assay Kits		
AF Cholesterol Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK436
AF HDL and LDL/VLDL Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK331
CETP Activity Assay Kit	sufficient for 100 colorimetric tests	» MAK106
CETP RP Activity Assay Kit	sufficient for 100 fluorometric tests	» MAK305
HMG-CoA Reductase Assay Kit	sufficient for 30 tests	» CS1090
LCFC-LCAT Activity Assay Kit	sufficient for 100 fluorometric tests	» MAK306
Lipase Assay Kit	sufficient for 100 colorimetric tests	» MAK482
LPL Activity Assay Kit	sufficient for 100 fluorometric tests	» MAK109
MTP Activity Assay Kit	sufficient for 100 fluorometric tests	» MAK110
PLTP Activity Assay Kit	sufficient for 100 fluorometric tests	» MAK108
Amino Acid Metabolism Assay Kits		
Alanine Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK500
ALT Activity Assay	sufficient for 100 colorimetric or fluorometric tests	» MAK052
Ammonia Assay Kit	sufficient for 100 tests	» AA0100
Ammonia Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK538
Arginase Activity Assay Kit	sufficient for 200 colorimetric tests	» MAK112
Arginase Assay Kit	sufficient for 100 colorimetric tests	» MAK533
Aspartate Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK495
Aspartate Transaminase (AST) Assay Kit	sufficient for 100 colorimetric tests	» MAK467

Name	Description	Cat No.
Branched Chain Amino Acid Kit	sufficient for 100 colorimetric tests	» MAK562
Creatine Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK494
Creatinine Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK475
D-Amino Acid Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK537
Formaldehyde Assay Kit	sufficient for 100 fluorometric tests	» MAK131
Glutamate Assay Kit	sufficient for 100 colorimetric tests	» MAK330
Glutamate Dehydrogenase (GDH) Assay Kit	sufficient for 100 colorimetric tests	» MAK499
Hydroxyproline Assay Kit	sufficient for 100 colorimetric tests, perchlorate-free	» MAK463
Indican Assay Kit	sufficient for 100 colorimetric tests	» MAK128
L-Amino Acid Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK490
Phenylalanine Assay Kit	sufficient for 100 fluorometric tests	» MAK484
Tryptophan Assay Kit	sufficient for 100 fluorometric tests	» MAK523
Urea Assay Kit III	sufficient for 100 colorimetric tests	» MAK471
Uric Acid Assay Kit	sufficient for 250 colorimetric tests	» MAK483
Xanthine Oxidase Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK497
Carbohydrate Metabolism Assay Kits		
Acetate Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK474
Mannitol Assay Kit	sufficient for 100 colorimetric tests	» MAK514
Sorbitol Assay Kit	sufficient for 100 colorimetric tests	» MAK442
Fructose Assay Kit	sufficient for 20 tests	» FA20
Glucose-6-Phosphate Dehydrogenase Assay Kit	sufficient for 100 colorimetric tests	» MAK451
Galactose Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK502
Glucose Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK476
Glucose (GO) Assay Kit	sufficient for 20 tests	» GAGO20
Glucose (HK) Assay Kit	sufficient for 20 tests	» GAHK20
Glucose Uptake Assay Kit	sufficient for 100 colorimetric tests	» MAK542
Glucose Uptake Assay Kit	sufficient for 100 fluorometric tests	» MAK489
Glucose-6-Phosphate Assay Kit	sufficient for 100 colorimetric tests	» MAK503
Glycogen Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK465

Name	Description	Cat No.
Glucose-6-Phosphate (G6P) Assay Kit	sufficient for 200 fluorometric tests	» MAK548
Lactose Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK487
Maltose Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK513
Pyruvate Dehydrogenase Activity Assay Kit	sufficient for 100 colorimetric tests	» MAK567
Sialic Acid Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK314
Sorbitol Dehydrogenase Assay Kit	sufficient for 100 colorimetric tests	» MAK317
Total Carbohydrate Assay Kit	sufficient for 100 colorimetric tests	» MAK559
α -Ketoglutarate Quantitation Kit	sufficient for 200 colorimetric tests	» MAK541
α -Amylase Activity Assay Kit	sufficient for 100 colorimetric tests	» MAK478
α -Glucosidase Activity Assay Kit	sufficient for 100 colorimetric tests	» MAK123
α -Mannosidase Assay Kit	sufficient for 100 colorimetric tests	» MAK318
β -Glucosidase Activity Assay Kit	sufficient for 100 colorimetric tests	» MAK129
Coenzymes and Cofactors Metabolism Assay Kits		
Acetyl-Coenzyme A Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK566
ADP Assay Kit	sufficient for 100 luminescent tests	» MAK133
ADP Assay Kit	sufficient for 100 fluorometric tests	» MAK518
ADP/ATP Ratio Assay Kit	sufficient for 100 luminescent tests	» MAK135
Coenzyme A Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK504
Fluorimetric cADP-Ribose Assay Kit	sufficient for 100 fluorometric tests	» MAK553
NAD/NADH Assay Kit	sufficient for 100 colorimetric tests	» MAK468
NADP/NADPH Assay Kit	sufficient for 100 colorimetric tests	» MAK479
Fatty Acid and Lipid Metabolism Assay Kits		
Acetate Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK474
Acetyl-Coenzyme A Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK566
Adipolysis Assay Kit	sufficient for 200 colorimetric or fluorometric tests	» MAK313
Fatty Acid Uptake Assay Kit	sufficient for 100 fluorometric tests	» MAK156
CETP Activity Assay Kit	sufficient for 100 fluorometric tests	» MAK106
Cholesterol Extraction Kit	sufficient for 40 extractions	» MAK175

Name	Description	Cat No.
Choline Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK508
Coenzyme A Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK504
Fatty Acid Extraction Kit	sufficient for 40 extractions	» MAK174
Fatty Acid Methylation Kit	sufficient for 100 tests	» MAK224
Fatty Acid Methyl Ester Purification Kit	sufficient for 50 tests	» MAK225
Free Fatty Acid Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK466
Free Glycerol Determination Kit	sufficient for 1000 tests	» FG0100
Glycerol Assay Kit	sufficient for 200 colorimetric or fluorometric tests	» MAK117
Adipolysis Assay Kit	sufficient for 200 colorimetric or fluorometric tests	» MAK313
Ketone Body Assay Kit	sufficient for 200 tests (UV)	» MAK134
Lipase Assay Kit	sufficient for 100 colorimetric tests	» MAK482
Lipid (Oil Red O) Staining Kit	sufficient for 2 96-well plates	» MAK560
Lipid Peroxidation (MDA) Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK568
LPL Activity Assay Kit	sufficient for 100 fluorometric tests	» MAK109
MTP Activity Assay Kit	sufficient for 100 fluorometric tests	» MAK110
Phospholipase D Activity Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK137
Phospholipid Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK122
High Sensitivity Triglyceride Fluorometric Assay Kit	sufficient for 100 fluorometric tests	» MAK565
High Sensitivity Beta-Hydroxybutyrate (Ketone Body) Assay Kit	sufficient for 200 fluorometric tests	» MAK547
PLTP Activity Assay Kit	sufficient for 100 fluorometric tests	» MAK108
Serum Triglyceride Determination Kit	sufficient for 250 tests	» TR0100
Triglyceride Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK564
Beta-Hydroxybutyrate (Ketone Body) Assay Kit	sufficient for 200 colorimetric tests	» MAK540
Inorganic Ions Metabolism Assay Kit		
Calcium Assay Kit	sufficient for 500 colorimetric tests	» MAK477
Fluo-4 No Wash Calcium Assay Kit	sufficient for 10 fluorometric test plates	» MAK552
Chloride Assay Kit	sufficient for 250 colorimetric tests	» MAK510
Copper Assay Kit	sufficient for 250 colorimetric tests	» MAK127

Name	Description	Cat No.
Iron Assay Kit	sufficient for 250 colorimetric tests	» MAK472
Magnesium Assay Kit	sufficient for 250 colorimetric tests	» MAK507
Nitrite/Nitrate Assay Kit, colorimetric	sufficient for 100 colorimetric tests	» 23479
Phosphate Assay Kit	sufficient for 500 colorimetric tests	» MAK488
Sulfate Assay Kits	sufficient for 200 turbidimetric tests	» MAK132
Zinc Assay Kit	sufficient for 250 colorimetric tests	» MAK496
Nutritional Analysis and Quantification Kits		
Adipolysis Assay Kit	sufficient for 200 colorimetric or fluorometric tests	» MAK313
Alcohol Dehydrogenase (ADH) Assay Kit	sufficient for 100 colorimetric tests	» MAK498
α -Amylase Activity Assay Kit	sufficient for 100 colorimetric tests	» MAK478
Antioxidant Assay Kit	sufficient for 100 colorimetric tests	» MAK334
ATP Assay Kit	sufficient for 100 assays (bioluminescent)	» MAK473
Autophagy Assay Kit	sufficient for 200 fluorometric tests	» MAK138
High Sensitivity Beta-Hydroxybutyrate (Ketone Body) Assay Kit	sufficient for 200 fluorometric tests	» MAK547
Beta-Lactamase Activity Assay Kit	sufficient for 200 colorimetric tests	» MAK545
β -Secretase (BACE1) Activity Detection Kit	sufficient for 250 fluorometric tests	» CS0010
Catalase Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK531
Citrate Synthase Assay Kit	sufficient for 480 colorimetric reactions	» CS0720
Colorimetric Aldehyde Assay Kit	sufficient for 200 colorimetric tests	» MAK139
Colorimetric Aldehyde Assay Kit, Blue	sufficient for 200 colorimetric tests	» MAK140
Colorimetric Biotin Assay Kit	sufficient for 200 colorimetric tests	» MAK171
Colorimetric Maleimide Assay Kit	sufficient for 100 colorimetric tests	» MAK162
Colorimetric Sphingomyelinase Assay Kit	sufficient for 200 colorimetric tests	» MAK152
Enterokinase (Enteropeptidase) Activity Assay Kit	sufficient for 200 colorimetric tests	» MAK544
Ethanol Assay Kit	sufficient for 500 colorimetric tests	» MAK480
Ethanol Assay Kit	sufficient for 100 colorimetric tests	» MAK481
Fatty Acid Extraction Kit	sufficient for 40 extractions	» MAK174
Fatty Acid Uptake Assay Kit	sufficient for 100 fluorometric tests	» MAK156
Ferric Reducing Antioxidant Power (FRAP) Assay Kit	sufficient for 250 colorimetric tests	» MAK509

Name	Description	Cat No.
Fluorimetric Hydrogen Peroxide Assay Kit	sufficient for 500 fluorometric tests (red fluorescence)	» MAK165
Fluorimetric Hydrogen Peroxide Assay Kit	sufficient for 500 fluorometric tests (infrared fluorescence)	» MAK166
Fluorometric Intracellular pH Assay Kit	sufficient for 1000 fluorometric tests	» MAK150
Fluorometric Intracellular ROS KIT	sufficient for 200 fluorometric tests (deep red)	» MAK142
Fluorometric Intracellular ROS KIT	sufficient for 200 fluorometric tests (green)	» MAK143
Fluorometric Intracellular ROS KIT	sufficient for 200 fluorometric tests (orange)	» MAK144
Fluorometric Intracellular ROS KIT	sufficient for 200 fluorometric tests (red)	» MAK145
Fluorometric Maleimide Assay Kit	sufficient for 200 fluorometric tests	» MAK167
Fluorometric Sphingomyelin Assay Kit	sufficient for 100 fluorometric tests	» MAK154
Fluorometric Thiol Assay Kit	sufficient for 200 fluorometric tests	» MAK151
Formate Assay Kit	sufficient for 100 colorimetric tests	» MAK491
Fructose Assay Kit	sufficient for 20 tests	» FA20
Fumarase Assay Kit	sufficient for 100 colorimetric tests	» MAK521
Galactose Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK502
Glucose Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK476
Glucose (GO) Assay Kit	sufficient for 20 tests	» GAGO20
Glucose (HK) Assay Kit	sufficient for 20 tests	» GAHK20
Glycerol 3-Phosphate (G3P) Assay Kit	sufficient for 200 colorimetric tests	» MAK546
Glycogen Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK465
High Sensitivity Glucose Quantitation Kit	sufficient for 500 fluorometric tests	» MAK543
Intracellular Hydrogen Peroxide Assay	sufficient for 200 fluorometric tests (green fluorescence)	» MAK164
Lactose Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK487
Lipid (Oil Red O) Staining Kit	sufficient for 2 96-well plates	» MAK560
Mannitol Assay Kit	sufficient for 100 colorimetric tests	» MAK514
Malate Dehydrogenase Assay Kit	sufficient for 100 colorimetric tests	» MAK512
Maltose Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK513
Mitochondrial Membrane Potential Kit	sufficient for 500 fluorometric tests (microplate readers)	» MAK159

Name	Description	Cat No.
Mitochondrial Membrane Potential Kit	sufficient for 100 fluorometric tests (flow cytometry)	» MAK160
Multidrug Resistance Assay Kit	Fluorometric MDR Assay	» MAK161
Oxalate Assay Kit	sufficient for 100 colorimetric tests	» MAK315
Proteasome 20S Activity Assay Kit	sufficient for 100 assays in 96 well plates	» MAK172
Pyrophosphate Assay Kit	sufficient for 200 fluorometric tests (Blue fluorescence)	» MAK168
High Sensitivity Pyrophosphate Assay Kit	sufficient for 200 fluorometric tests	» MAK169
Pyruvate Dehydrogenase Activity Assay Kit	sufficient for 100 colorimetric tests	» MAK567
Renin Assay Kit	sufficient for 100 fluorometric tests	» MAK157
Sorbitol Assay Kit	sufficient for 100 colorimetric tests	» MAK442
Starch Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK522
Starch (GO/P) Assay Kit	sufficient for 20 tests	» STA20
Starch (HK) Assay Kit	sufficient for 20 tests	» SA20
Succinate Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK335
Succinate Dehydrogenase Assay Kit	sufficient for 100 colorimetric tests	» MAK561
Sucrose Assay Kit	sufficient for 20 tests	» SCA20
Total Dietary Fiber Assay Kit	sufficient for 100 tests	» TDF100A
Total Dietary Fiber Assay Control Kit	sufficient for 10 assays	» TDFC10
Triglyceride Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK564
High Sensitivity Triglyceride Fluorometric Assay Kit	sufficient for 100 fluorometric tests	» MAK565
Tryptophan Assay Kit	sufficient for 100 fluorometric tests	» MAK523
Universal Fluorometric Kinase Assay Kit	sufficient for 250 fluorometric tests	» MAK173
Blood and Urine Chemistry Assay Kits		
Acetate Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK474
Ketone Body Assay	sufficient for 200 tests (UV)	» MAK134
Acid Phosphatase Fluorometric Assay Kit	sufficient for 100 fluorometric tests	» MAK525
Alanine Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK500
Alcohol Dehydrogenase (ADH) Assay Kit	sufficient for 100 colorimetric tests	» MAK498
ALT Activity Assay	sufficient for 100 colorimetric or fluorometric tests	» MAK052
Ammonia Assay Kit	sufficient for 100 tests	» AA0100
Ammonia Assay Kit	sufficient for 200 fluorometric tests	» MAK310

Name	Description	Cat No.
Ammonia Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK538
α-Amylase Activity Assay Kit	sufficient for 100 colorimetric tests	» MAK478
Arginase Activity Assay Kit	sufficient for 200 colorimetric tests	» MAK112
Arginase Assay Kit	sufficient for 100 colorimetric tests	» MAK533
Ascorbic Acid Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK505
Aspartate Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK495
Aspartate Transaminase (AST) Assay Kit	sufficient for 100 colorimetric tests	» MAK467
BCG Albumin Assay Kit	sufficient for 250 colorimetric tests	» MAK124
BCP Albumin Assay Kit	sufficient for 250 colorimetric tests	» MAK125
Beta-Hydroxybutyrate (Ketone Body) Assay Kit	sufficient for 200 colorimetric tests	» MAK540
Bile Acid Assay Kit	sufficient for 100 fluorometric tests	» MAK309
Bilirubin Assay Kit	sufficient for 180 colorimetric tests	» MAK126
Butyrylcholinesterase (BChE) Activity Assay Kit	sufficient for 100 colorimetric tests	» MAK551
Calcium Assay Kit	sufficient for 500 colorimetric tests	» MAK477
Chloride Assay Kit	sufficient for 250 colorimetric tests	» MAK510
AF Cholesterol Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK436
Choline Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK508
Citrate Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK333
Coenzyme A Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK504
Copper Assay Kit	sufficient for 250 colorimetric tests	» MAK127
Creatine Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK494
Creatine Kinase Activity Assay Kit	sufficient for 100 UV tests	» MAK116
Creatinine Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK475
Ethanol Assay Kit	sufficient for 500 colorimetric tests	» MAK480
Ethanol Assay Kit	sufficient for 100 colorimetric tests	» MAK481
Formaldehyde Assay Kit	sufficient for 100 fluorometric tests	» MAK131
Free Fatty Acid Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK466

Name	Description	Cat No.
Free Glycerol Determination Kit	sufficient for 1000 tests	» FG0100
Glutamate Assay Kit	sufficient for 100 colorimetric tests	» MAK330
Glutamate Dehydrogenase (GDH) Assay Kit	sufficient for 100 colorimetric tests	» MAK499
Glutathione (GSH) Assay Kit	sufficient for 250 colorimetric tests	» MAK517
Glutathione Reductase Kit	sufficient for 100 colorimetric tests	» MAK535
Glycerol Assay Kit	sufficient for 200 colorimetric or fluorometric tests	» MAK117
AF HDL and LDL/VLDL Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK331
Heme Assay Kit	sufficient for 250 colorimetric tests	» MAK316
Hemoglobin Assay Kit	sufficient for 250 colorimetric tests	» MAK115
Indican Assay Kit	sufficient for 100 colorimetric tests	» MAK128
Iron Assay Kit	sufficient for 250 colorimetric tests	» MAK472
Ketone Body Assay Kit	sufficient for 200 UV tests	» MAK134
L-Lactate Assay Kit	sufficient for 100 colorimetric tests	» MAK329
L-Lactate Assay Kit	sufficient for 100 fluorometric tests	» MAK443
Lactate Dehydrogenase Assay Kit	sufficient for 100 colorimetric tests	» MAK464
Lipase Assay Kit	sufficient for 100 colorimetric tests	» MAK482
Magnesium Assay Kit	sufficient for 250 colorimetric tests	» MAK507
Malachite Green Phosphate Assay Kit	sufficient for 2500 colorimetric tests	» MAK307
Monoamine Oxidase Activity Assay Kit	sufficient for 100 fluorometric tests	» MAK136
Oxalate Assay Kit	sufficient for 100 colorimetric tests	» MAK315
Oxaloacetate Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK515
Peroxide Assay Kit	sufficient for 250 colorimetric tests	» MAK311
Phenylalanine Assay Kit	sufficient for 100 fluorometric tests	» MAK484
Phosphate Assay Kit	sufficient for colorimetric tests	» MAK308
Phosphate Assay Kit	sufficient for 500 colorimetric tests	» MAK488
Phospholipid Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK122
Pyruvate Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK332
Salicylate Assay Kit	sufficient for 100 colorimetric tests	» MAK534
Serum Triglyceride Determination Kit	sufficient for 250 tests	» TR0100

Name	Description	Cat No.
Sulfate Assay Kits	sufficient for 200 turbidimetric tests	» MAK132
Tryptophan Assay Kit	sufficient for 100 fluorometric tests	» MAK523
Urea Assay Kit III	sufficient for 100 colorimetric tests	» MAK471
Zinc Assay Kit	sufficient for 250 colorimetric tests	» MAK496
Enzyme Activity Assay Kits		
Acetylcholinesterase Activity Assay Kit	sufficient for 100 colorimetric tests	» MAK119
Fluorimetric Acetylcholinesterase Assay Kit	sufficient for 200 fluorometric tests	» MAK554
Acid Phosphatase Fluorometric Assay Kit	sufficient for 100 fluorometric tests	» MAK525
Aconitase Assay Kit	sufficient for 100 colorimetric tests	» MAK337
Alcohol Dehydrogenase (ADH) Assay Kit	sufficient for 100 colorimetric tests	» MAK498
Alkaline Phosphatase (ALP) Assay Kit	sufficient for 100 fluorometric tests	» MAK530
ALT Activity Assay	sufficient for 100 colorimetric or fluorometric tests	» MAK052
Arginase Activity Assay Kit	sufficient for 200 colorimetric tests	» MAK112
Arginase Assay Kit	sufficient for 100 colorimetric tests	» MAK533
Aspartate Transaminase (AST) Assay Kit	sufficient for 100 colorimetric tests	» MAK467
ATPase/GTPase Activity Assay Kit	sufficient for 200 colorimetric tests	» MAK113
Beta-Lactamase Activity Assay Kit	sufficient for 200 colorimetric tests	» MAK545
β -Secretase (BACE1) Activity Detection Kit	sufficient for 250 fluorometric tests	» CS0010
Butyrylcholinesterase (BChE) Activity Assay Kit	sufficient for 100 colorimetric tests	» MAK551
Catalase Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK531
Diamine Oxidase Assay Kit	sufficient for 100 fluorometric tests	» MAK526
Enterokinase (Enteropeptidase) Activity Assay Kit	sufficient for 200 colorimetric tests	» MAK544
Glucose Oxidase Activity Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK501
α -Glucosidase Activity Assay Kit	sufficient for 100 colorimetric tests	» MAK123
β -Glucosidase Activity Assay Kit	sufficient for 100 colorimetric tests	» MAK129
Glutamate Dehydrogenase (GDH) Assay Kit	sufficient for 100 colorimetric tests	» MAK499
Glyoxalase I Activity Assay Kit	sufficient for 100 colorimetric tests	» MAK114
Isocitrate Dehydrogenase (IDH) Assay Kit	sufficient for 100 colorimetric tests	» MAK493
Invertase Activity Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK118

Name	Description	Cat No.
Lactate Dehydrogenase Assay Kit	sufficient for 100 colorimetric tests	» MAK464
LDH Cytotoxicity Assay Kit	sufficient for 100 colorimetric tests	» MAK529
Lipase Assay Kit	sufficient for 100 colorimetric tests	» MAK482
LPL Activity Assay Kit	sufficient for 100 fluorometric tests	» MAK109
Lysyl Oxidase (LOX) Assay Kit	sufficient for 500 fluorometric tests	» MAK555
Monoamine Oxidase Activity Assay Kit	sufficient for 100 fluorometric tests	» MAK136
Monoamine Oxidase (MAO) Inhibitor Screening Kit	sufficient for 100 fluorometric tests	» MAK520
MPO Colorimetric Activity Assay Kit	sufficient for 100 colorimetric tests	» MAK563
MPO Fluorometric Activity Assay Kit	sufficient for 100 fluorometric tests	» MAK516
MTP Activity Assay Kit	sufficient for 100 fluorometric tests	» MAK110
Neuraminidase Activity Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK121
Nitric Oxide Synthase (NOS) Assay Kit	sufficient for 100 colorimetric tests	» MAK532
Peroxidase Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK506
Phospholipase D Activity Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK137
PLTP Activity Assay Kit	sufficient for 100 fluorometric tests	» MAK108
Pyruvate Kinase Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK485
Succinate Dehydrogenase Assay Kit	sufficient for 100 colorimetric tests	» MAK561
Superoxide Dismutase (SOD) Assay Kit	sufficient for 100 colorimetric tests	» MAK528
Trypsin Activity Assay Kit	sufficient for 100 colorimetric tests	» MAK558
Tyrosinase Assay Kit	sufficient for 100 colorimetric tests	» MAK550
Urease Activity Assay Kit	sufficient for 100 colorimetric tests	» MAK120
Xanthine Oxidase Assay Kit	sufficient for 100 colorimetric or fluorometric tests	» MAK497
Extraction Kits		
Fatty Acid Extraction Kit	sufficient for 40 extractions	» MAK174
Cholesterol Extraction Kit	sufficient for 40 extractions	» MAK175

To place an order or receive technical assistance

Order/Customer Service:
SigmaAldrich.com/order

Technical Service:
SigmaAldrich.com/techservice

Safety-related Information:
SigmaAldrich.com/safetycenter

SigmaAldrich.com

For more information,
visit **SigmaAldrich.com/assaykits**

Merck KGaA
Frankfurter Strasse 250
64293 Darmstadt, Germany

We have built a unique collection of life science brands with
unrivalled experience in supporting your scientific advancements.

Millipore® Sigma-Aldrich® Supelco® Milli-Q® SAFC® BioReliance®

