

Product Information

Conductive Single-Wall Carbon Nanotubes

Catalog Number **704113**
 Store at Room Temperature

Synonym: SWCNT

Product Description

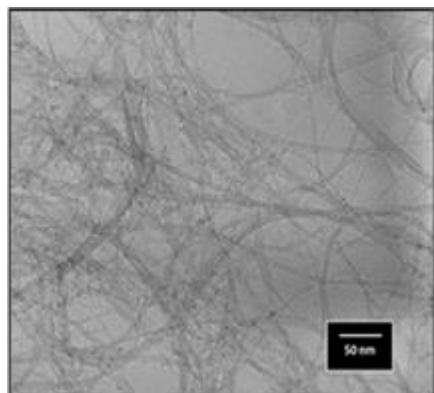
These Single-Wall Carbon Nanotubes (SWCNT) are produced using patented CoMoCAT™ synthesis technology. 704113 is the most economical SWCNT product, offering good electrical conductivity and high carbon purity at an affordable price.

Representative product characteristics are presented.

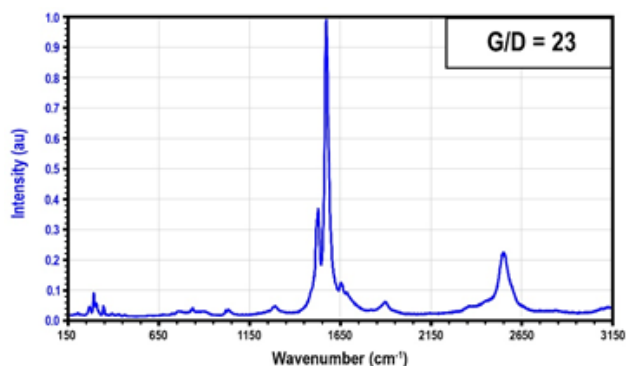
Typical Product Properties

Property (Measurement)	704113
Carbon Purity (TGA)	≥90 wt%
CNT Purity (TGA)	≥89 wt%
Average Diameter (NIRF)	0.82 nm
Median Length (AFM)	1 μm
Bulk Density (ASTM D7481)	0.1 g/cm ³
Moisture Content (TGA)	≤5 wt%
Specific Surface Area (BET)	≥700 m ² /g
G/D Ratio (Raman [633 nm])	≥15
Standard Product Form	Powder

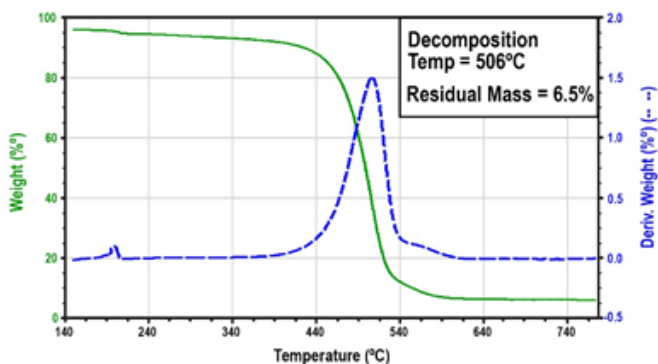
Typical SEM



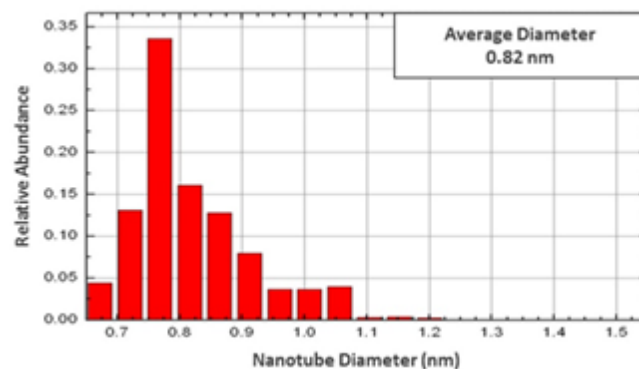
Typical Raman Analysis (633 nm)



Typical Thermogravimetric Analysis

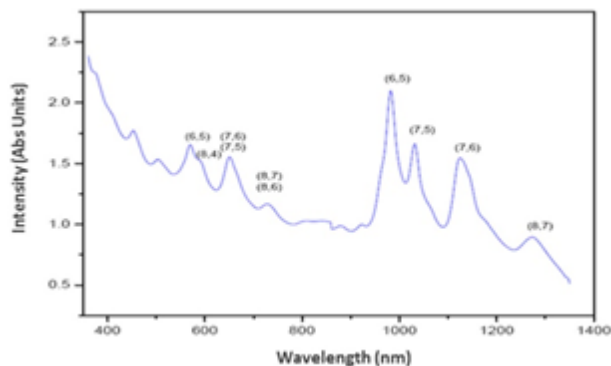


Typical Diameter Distribution



Typical Optical Absorbance

OA Peaks Corresponding to Eii Transition Energies of Included Chiral Species



Precautions and Disclaimer

This product is for R&D use only, not for drug, household, or other uses. Please consult the Safety Data Sheet for information regarding hazards and safe handling practices.

CoMoCAT is a trademark of Chasm Advanced Materials, Inc.

RC,MAM 05/17-1