

## Product Information

### (6,5) Enriched Single-Wall Carbon Nanotubes

Catalog Number **773735**  
 Store at Room Temperature

Synonym: SWCNT

#### Product Description

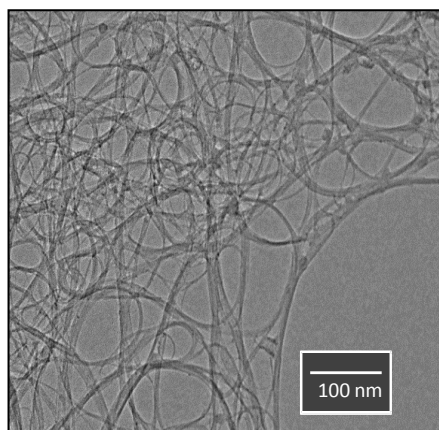
These Single-Wall Carbon Nanotubes (SWCNT) are produced using patented CoMoCAT™ synthesis technology. 773735 is a single-wall carbon nanotube product uniquely enriched in semiconducting tubes, with (6,5) chirality being the most abundant. ~95% of the SWCNTs in 773735 are semiconducting, with ~41% of those tubes being (6,5) chirality. 773735 is targeted for transistor, sensor, and biomedical applications.

Representative product characteristics are presented.

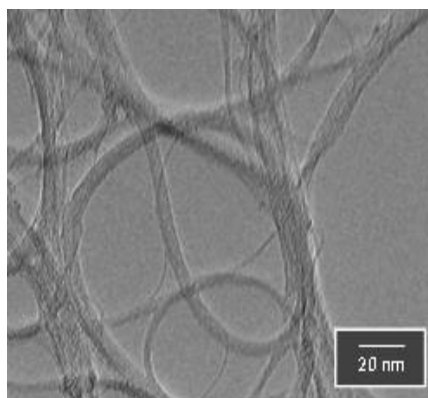
#### Typical Product Properties

Property (Measurement)	773735
Carbon Purity (TGA)	≥95 wt%
CNT Purity (TGA)	≥94 wt%
Semiconducting CNT Content (Opt. Abs)	95%
(6,5) CNT Content (NIRF)	≥40%
Average Diameter (NIRF)	0.78 nm
Median Length (AFM)	1 μm
Bulk Density (ASTM D7481)	0.1 g/cm <sup>3</sup>
Moisture Content (TGA)	≤5 wt%
Specific Surface Area (BET)	≥700 m <sup>2</sup> /g
G/D Ratio (Raman [633 nm])	≥20
Standard Product Form	Powder

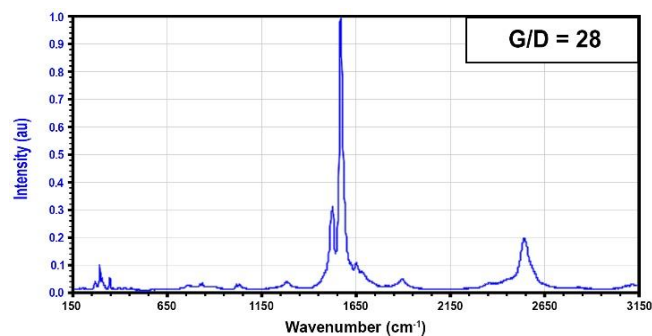
#### Typical SEM



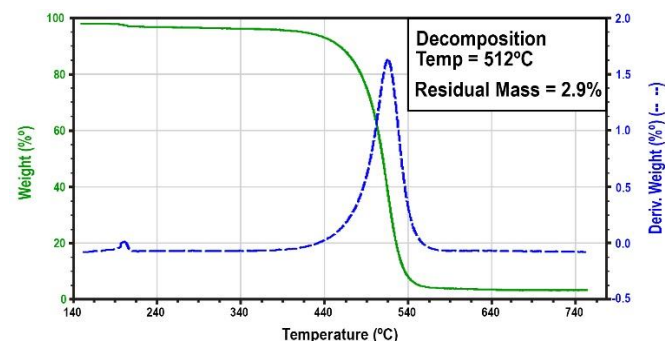
#### Typical SEM



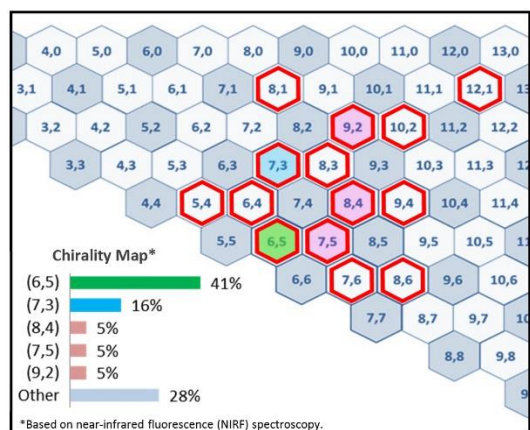
#### Typical Raman Analysis (633 nm)



#### Typical Thermogravimetric Analysis

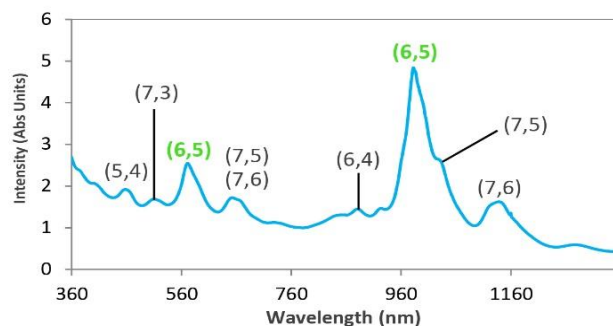


### Typical Chirality Distribution

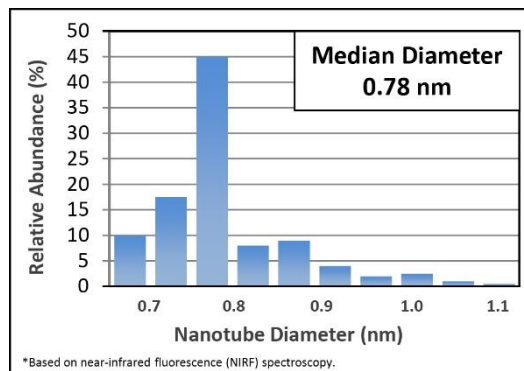


### Typical Optical Absorbance

OA Peaks Corresponding to Eii Transition Energies of Included Chiral Species



### Diameter Distribution



### 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