

## Product Information

### Large Diameter Single-Wall Carbon Nanotubes

Catalog Number **724777**  
 Store at Room Temperature

Synonym: SWCNT

#### Product Description

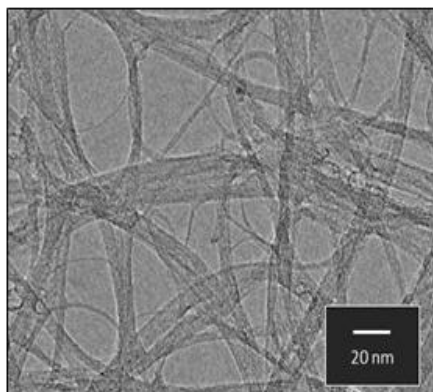
These Single-Wall Carbon Nanotubes (SWCNT) are produced using patented CoMoCAT™ synthesis technology. 724777 is the largest-diameter SW CNT product.

Representative product characteristics are presented.

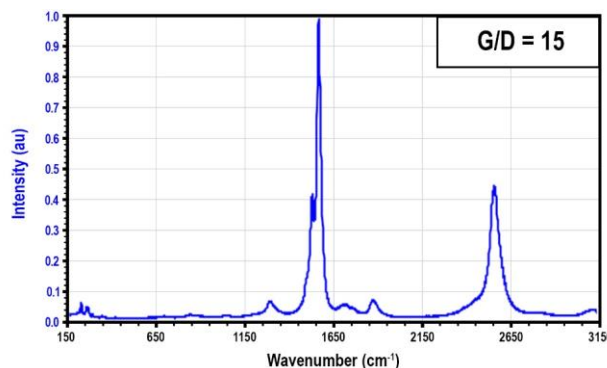
#### Typical Product Properties

Property (Measurement)	724777
Carbon Purity (TGA)	≥90 wt%
CNT Purity (TGA)	≥89 wt%
Average Diameter (NIRF)	1.3 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])	≥15
Standard Product Form	Powder

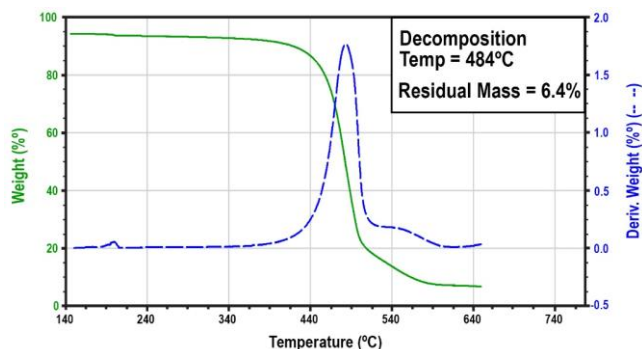
#### Typical SEM



#### Typical Raman Analysis (633 nm)



#### Typical Thermogravimetric Analysis



#### 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