

big flask even bigger results

The Thermo Scientific™ Nunc™ 300 cm² Cell Culture Flask is our newest and biggest cell culture flask. In fact, this 1,900 mL total volume flask provides the largest culture area of any single-layer flask on the market, and is cell culture treated to deliver a uniform surface chemistry for promoting cell attachment and expanded growth. It's another example of the Thermo Fisher Scientific commitment to advancing cell culture research.



high-volume cell growth



The Nunc™ 300 cm² flask expands upon our previous product range of 25 cm² - 225 cm² flasks, dishes and multidishes.



The new **Thermo Scientific Nunc 300 cm² Cell Culture Flask** offers cell culture researchers numerous workflow benefits.

- Promotes cell attachment and growth**
 Vacuum-gas plasma tissue-culture treated polystyrene for uniform surface chemistry, cell attachment, spreading and promoting cell growth
- Durable construction, friendly design**
 Blow-molded, one-piece design with straight neck and molded grip notches that make this 1,900 mL flask easy to hold
- Quick volume estimation and sample identification**
 Includes bold printed graduations – from 100 mL to 500 mL, in 50 mL graduations – and large writing patch
- Maintains optimal sterility during growth and passage**
 Filter cap incorporates a 0.2 µm pore size hydrophobic membrane that allows gas exchange while maintaining sterility
- High-quality assurance**
 Sterilized by gamma irradiation, non-pyrogenic
- Easy to stack**
 Prominent stacking feet on upper surface facilitate reliable stacking in incubators and under your culture hood

Ordering Information

Catalog number	Surface	Closure	Culture Area, cm ²	Working Volume, mL	No. per Pack	No. per Case
132097	Treated	Vent/close	300	1,900	1	12
132098	Treated	Filter	300	1,900	1	12

© 2012 Thermo Fisher Scientific, Inc. All rights reserved. All other trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries.

ANZ: Australia: 1300 735 292, New Zealand: 0800 933 966 **Asia:** China Toll-free: 800-810-5118 or 400-650-5118; India: +91 22 6716 2200, India Toll-free: 1 800 22 8374; Japan: +81 3 5826 1616; Other Asian countries: 65 68729717
Europe: Austria: +43 1 801 40 0; Belgium: +32 53 73 42 41; Denmark: +45 4631 2000; France: +33 2 2803 2180; Germany: +49 6184 90 6940, Germany Toll-free: 08001-536 376; Italy: +39 02 02 95059 or 434-254-375; Netherlands: +31 76 571 4440; Nordic/Baltic countries: +358 9 329 100; Russia/CIS: +7 (812) 703 42 15; Spain/Portugal: +34 93 223 09 18; Switzerland: +41 44 454 12 12; UK/Ireland: +44 870 609 920
North America: USA/Canada +1 585 586 8800; USA Toll-free: 800 625 4327 **South America:** USA sales support: +1 585 899 7198
Countries not listed: +49 6184 90 6940 or +33 2 2803 2180

www.thermoscientific.com/cellcultureflasks

Thermo
SCIENTIFIC

Part of Thermo Fisher Scientific