

FNC Coating Mix®

FNC Coating Mix® is formulated to dramatically increase the attachment of cells to the plastic growth substratum used for culturing mammalian cell lines. Many fastidious cell types such as human prostate and breast epithelial cells plated in the appropriate serum-free medium will attach and grow more effectively on FNC-coated plastic surfaces.

Ordering Information

[Order Online »](#)

AES-0407-100	100ml
--------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability 6 months at +4°C

Handling Do not freeze.

Short Term Storage +4°C

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Formulation Liquid; Ready-to-Use.

Technical Info / Product Notes **pH:** 7.1 -7.4
Osmolality: 280 -300mOsM

For the Original Manufacturer's data sheet please [click here](#).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com