

SILAC – DMEM/F12

DMEM/F12 is a serum-free medium formulation for general use. This formulation is a 1:1 blend of DMEM and Ham's F12 media supplied complete, ready-to-use with L-glutamine, Hepes, BPE and EGF for culturing a wide range of cell types. Contains no phenol red or antibiotics.

Ordering Information

[Order Online »](#)

AES-0423	500ml
----------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability Stability at 4°C: 4-6 weeks.

Short Term Storage +4°C

Long Term Storage -80°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Formulation Liquid; Ready-to-Use.



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