

EGF receptor (human)

EGFR is active but not autophosphorylated.

Citations: 4

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-SE116-0010	10U
----------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Epidermal growth factor receptor, EGFR
Formulation	Liquid. In 10% glycerol, 0.05% Triton X-100 buffer containing 5 mM EGF.
MW	170 kDa
Purity	≥90% (SDS-PAGE)
Source	Isolated from human A431 cells.
Specific Activity	One unit will add one pmol phosphate to angiotensin II per minute at pH 7.4 and 30°C
UniProt ID	P00533

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

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

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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