

c-erbB-2 (human), (recombinant)

c-erbB2 is best known in its role in breast cancer, where it is amplified in approximately 20% of invasive cancers. The humanized monoclonal antibody, trastuzumab, and the small molecule, lapatinib, inhibit c-erbB2.

Ordering Information

[Order Online »](#)

BML-SE460-0010	10µg
-----------------------	------

Manuals, SDS & CoFA

[View Online »](#)

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name ErbB2, Her2, Neu, CD340

Application Notes Useful for kinetic and functional studies, phosphorylation of target substrates, drug screening.

MW 92 kDa

Source Produced in insect cells. Active c-erbB2 (aa 679-1255).
Produced in a baculovirus expression system.

UniProt ID P04626

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)