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



eu@enzolifesciences.com