

14-3-3 γ (human), (recombinant) (GST- tag)

14-3-3 proteins are dimeric phosphoserine/phosphothreonine binding proteins, which have been shown to be involved in the regulation of many crucial processes such as signal transduction, cell-cycle control, apoptosis, protein localization, phosphorylation-state, stability, transcription, metabolism, and malignant transformation. Potential 14-3-3 γ protein–protein interactions are also differentially regulated by cytokine stimulation. One report showed that 14-3-3 γ expression is induced in arterial trauma by cytokines, which suggests that this protein may play an important role in progression of vascular proliferative diseases.

Citations: 2

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-SE313-0100	100 μ g
----------------	-------------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Notes Can be used for protein-protein interaction studies and gel overlays.

Formulation Liquid. In 50mM TRIS/HCl, pH 7.5, 150mM NaCl.

MW ~60 kDa

Purity ≥90% (SDS-PAGE)

Purity Detail Purified by multi-step chromatography.

Source Produced in *E. coli*. Human 14-3-3γ is fused at the N-terminus to a GST-tag.

UniProt ID P61981



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)