

# 14-3-3 $\epsilon$ (human), (recombinant) (GST- His-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 and stability, transcription, metabolism, and malignant transformation. 14-3-3 $\epsilon$  is implicated in an early left-right patterning in amphibian embryos. In addition, children with Miller-Dieker syndrome, a severe brain malformation associated with mental retardation, have heterozygous deletions of both the 14-3-3 $\epsilon$  and LIS1 genes. It has also been shown to be the only 14-3-3 isoform associated with prion protein amyloid deposits of Gerstmann-Straussler-Scheinker disease.

## Ordering Information

[Order Online »](#)

|                 |             |
|-----------------|-------------|
| ENZ-PRT208-0100 | 100 $\mu$ g |
|-----------------|-------------|

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

**Handling** Avoid freeze/thaw cycles.

**Long Term Storage** -80°C

**Shipping** Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Application Notes** Drug screening, protein-protein interaction studies, gel overlays.

**Formulation** Liquid. In 20 mM Tris-HCl, pH 8.0, containing 100 mM NaCl.

**MW** ~56 kDa

**Purity** ≥90% (SDS-PAGE)

**Purity Detail** Purified by multi-step chromatography.

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

**UniProt ID** P62258

**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)