

# 14-3-3 $\zeta$ (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 and stability, transcription, metabolism, and malignant transformation. 14-3-3  $\zeta$  has been shown to be present in brain amyloid plaques.

**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 <i>E. coli</i> . Human 14-3-3ζ is fused at the N-terminus to a GST-tag.
UniProt ID	P63104

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)