

14-3-3 ϵ (human), (recombinant) (untagged)

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 ϵ 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 ϵ 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.

Citations: 1

[View Online »](#)

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, containing 100 mM NaCl.
Purity	≥90% (SDS-PAGE)
Purity Detail	Purified by multi-step chromatography
Sequence	~30 kDa
Source	Produced in <i>E. coli</i> .
UniProt ID	P62258

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com