

14-3-3 θ /t (human), (recombinant) (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 θ /t has been shown to be present in testis and 14-3-3 θ mRNA is up-regulated in amyotrophic lateral sclerosis spinal cord. Both PKC μ and Bax have been shown to be negatively regulated by 14-3-3 θ

Ordering Information

[Order Online »](#)

BML-SE487-0100	100 μ g
----------------	-------------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -80°C.
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 50mM NaH ₂ PO ₄ , pH 8.0, 300mM NaCl.
MW	~30 kDa
Purity	≥90% (SDS-PAGE)
Purity Detail	Purified by multi-step chromatography.
Source	Produced in <i>E. coli</i> . Human 14-3-3 τ /θ is fused at the C-terminus to a His-tag.
UniProt ID	P27348



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)