

14-3-3 σ (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 σ is the main 14-3-3 isoform that has been directly implicated in the development of cancer. The 14-3-3 σ gene is directly regulated by p53 and is silenced by CpG methylation in a large number of carcinomas.

Manuals, SDS & CofA

[View Online »](#)

kDa

250

148

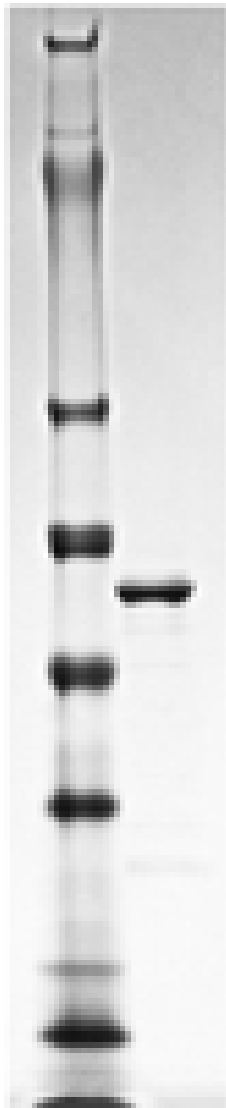
98

64

50

36

22



SDS-PAGE Analysis: Lane1: MWM; Lane 2: 1.0 μg of purified human 14-3-3 σ protein.

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 14-3-3σ is fused at the N-terminus to a GST-tag.
UniProt ID	P31947

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