

p62 (human), (recombinant) (GST- tag)

The p62 protein functions as a ubiquitin (Ub)-binding scaffold, which regulates a diverse range of signalling pathways leading to activation of the nuclear factor kappa B (NF-κB) family of transcription factors. p62 also plays an important role in the control of induced osteoclastogenesis, and mutations affecting the SQSTM1 gene are commonly found in patients with the skeletal disorder Paget's disease of bone (PDB). The vast majority of the PDB associated mutations identified cluster within the UBA domain, impairing p62's ability to bind Ub and resulting in dysregulated NF-κB signalling.

p62 is a critical regulator of the degradation of ubiquitinated proteins by macroautophagy. The p62 protein has a domain structure consistent with its participation in multiple signalling complexes, including a C-terminal Ub-associated (UBA) domain through which p62 binds non-covalently to Ub and a LC3 interacting region (LIR), facilitating simultaneous binding to ubiquitinated proteins and the autophagic machinery *via* LC3/GABARAP. As a result p62 acts as an important substrate receptor for selective autophagy.

Citations: 2

[View Online »](#)

Ordering Information

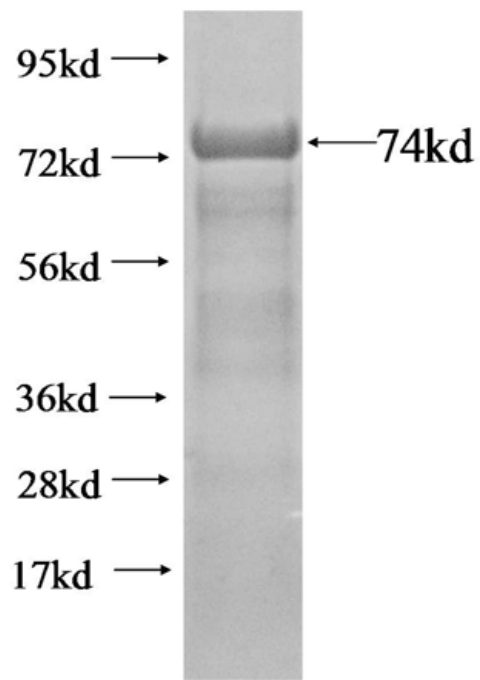
[Order Online »](#)

ENZ-PRT120-0050

50µg

Manuals, SDS & CofA

[View Online »](#)



Western Blot analysis of ENZ-PRT120.

Handling & Storage

Use/Stability	Stable for at least 12 months after receipt when stored at -20°C.
Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.
Short Term Storage	-20°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Sequestosome-1, EBI3-associated protein of 60 kDa
Formulation	Lyophilized from sterile PBS, pH 7.4, 5% trehalose, 5% mannitol.
MW	~74.4kDa
Purity	≥75% (SDS-PAGE)
Source	Produced in <i>E. coli</i> . Human p62 (1-440 aa) is fused to a N-terminal GST-tag.
UniProt ID	Q13501

Last modified: July 9, 2025



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)