

Proteasome activator 11S γ subunit (human), (recombinant)

A nuclear homo-heptameric form of the 11S proteasome activator built up by 11S subunit γ (designated 11S γ) was found to facilitate the degradation of an intact proteasome substrate by the 20S proteasome.

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-PW9875-0100	100 μ g
-----------------	-------------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Proteasome activator complex subunit 3, PA28γ

Formulation Liquid. In PBS pH 7.2 , 1mM DTT

MW ~29.5 kDa

Source Produced in *E.coli*.

UniProt ID P61289



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