

# Saposin C (human), (recombinant)

An essential activator for  
glucocerebrosidase, the enzyme deficient in  
Gaucher disease.

Glycoprotein with a variety of functions. Important for exchange of lipids  
from CD1 molecules to lysosomal antigens.

Citations: 4

[View Online »](#)

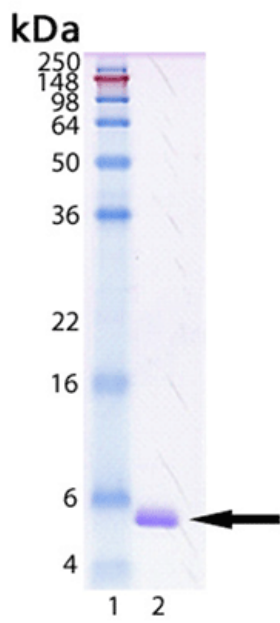
Ordering Information

[Order Online »](#)

ALX-201-262-C050	50µg
------------------	------

Manuals, SDS & CofA

[View Online »](#)



SDS-PAGE Analysis: Lane 1: MWM; Lane 2: 1.0  $\mu\text{g}$  of purified Saposin C.

## Handling & Storage

Use/Stability	Stable for at least one year after receipt when stored at +4°C.
Long Term Storage	+4°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	Co-β-glucosidase, A1 activator, Glucosylceramidase activator, Sphingolipid activator protein 2, SAP-2
Formulation	Lyophilized. In 50mM Tris, pH 6.0, containing 150mM sodium chloride.
MW	~8kDa
Purity	≥95% (SDS-PAGE)
Reconstitution	Reconstitute with 100µl sterile water.
Source	Produced in <i>E. coli</i> . Human recombinant saposin C (aa 309-388) is fused at the C-terminus to a His-tag and at the N-terminus to a T-7-tag.
UniProt ID	P07602

Last modified: May 29, 2024



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)