

Data Sheet

SIRP- α (CD172a), His-Tag

Human, recombinant, C-terminal His-tag

Catalog #: 71145

Lot #: 140331

Conc.: 2.15 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.4 mM KCl, and 20% glycerol.

Stability: >6 months at -80°C . Avoid freeze-thaw cycles. *Protein may be diluted to $\geq 100 \mu\text{g/ml}$ in PBS + glycerol and stored at -80°C .*

References:

1. Hatherley, D., *et al.*, *Mol. Cell*, 2008, **31**:266-77
2. Baba, N., *et al.*, *J. Exp. Med.*, 2013, **210**(6):1251-63

Description: Signal regulatory protein-alpha (SIRP- α), also known as Tyrosine-protein phosphatase non-receptor type substrate 1, CD172a, Bit, and MyD-1, GenBank Accession No. NM_080792, a.a. 27-370, with C-terminal His-tag, MW 38.4 kDa*, expressed in a HEK293 expression system. This protein acts as a receptor for CD47; a.a. 27-370 correspond to the extracellular portion of SIRP- α . *Note: SIRP- α is heavily glycosylated, resulting in apparent higher molecular weight.

Application: Useful for studying protein binding and screening small molecules.

Quality Assurance

4-20% SDS-PAGE **Coomassie staining**

Lane 1:

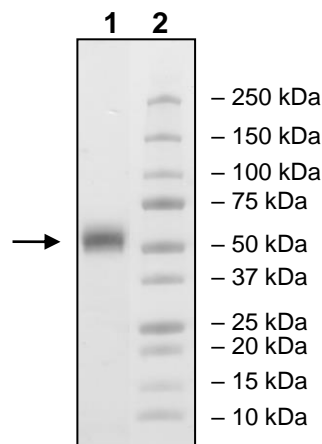
2.7 μg SIRP- α

Lane 2:

Protein Marker

MW: 38.4 kDa + glycans

Purity: $\geq 90\%$



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.259.5734**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6044 Cornerstone Court West, Suite E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Protein Sequence

GVAGEEELQVIQPDKSVLVAAGETATLRCTATSLIPVGPIQWFRGAGPGRELIYNQKEG
HFPRVTTVSDLTKRNNMDFSIRIGNITPADAGTYCYVKFRKGGSPDDVEFKSGAGTELSV
RAKPSAPVVSGPAARATPQHTVSFTCESHGFSPRDITLKWFKNGNELSDFQTNVDPVG
ESVSYSIHSTAKVVLTRDVDHSQVICEVAHVTLQGDPLRGTANLSETIRVPPTLEVTTQQP
VRAENQVNVTCQVRKFYPQRLQLTWLENGNVSRTETASTVTENKDGTYNWMSWLLVN
VSAHRDDVKLTCQVEHDGQPAVSKSHDLKVS AHPKEQGSNTAAENTGSNERHHHHHH

Black: SIRP α (a.a. 27-370)

Red: His(6x)

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.259.5734**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com