

Data Sheet

SMURF 1

Human, recombinant, N-terminal FLAG tag

Catalog #: 80402

Lot#: 120103

Conc.: 0.2 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0,
110 mM NaCl, 2.2 mM KCl, 80 ng/ul Flag
peptide, 3mM DTT, and 20% Glycerol.

Stability: >6 months at -80°C

References:

1. Narimatsu M. *et al.*, *Cell* **137** (2), 295-307 (2009)
2. Fukunaga E. *et al.*, *J. Biol. Chem.* **283** (51), 35660-35667 (2008).

Description:

Human SMURF 1 (GenBank Accession No. NM_020429), amino-acids 150-end, with N-terminal FLAG tag, MW=70 kDa, expressed in a Baculovirus infected Sf9 cell expression system.

Assay Conditions:

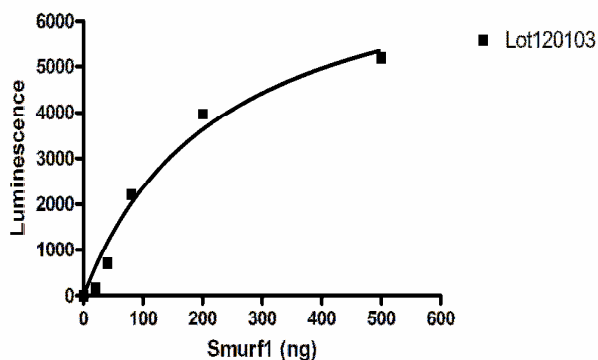
50 µl of reaction mix (4 mM ATP, 5 mM MgCl₂, 1 mM DTT, E1 (0.5 µg), E2 (2.8 µg), Biotin-Ubiquitin (0.5 µg) and Smurf 1 (0-1.25µg) in 50 mM Tris pH 7.4) was incubated for 30 min at 37°C. The reaction mixture was transferred to a streptavidin-coated plate and incubated for additional 30 min at 37°C. Anti-Smurf 1 and HRP-labeled secondary antibody were incubated with the reaction mixture for 1 hour each, followed by luminescent detection.

Application:

Useful in conjunction with E1 and E2 for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Assurance

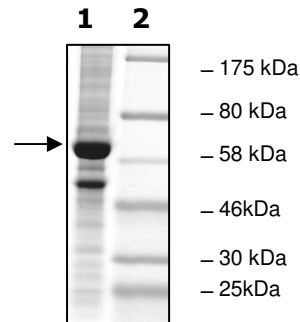
SMURF 1 Specific Activity



10% SDS-PAGE Coomassie staining

Lane 1:
3 µg SMURF 1
Lane 2:
Protein Marker
BioLabs (#P7708L)

MW: 70 kDa
Purity: ≥53%



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 **1.858.481.8694**

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

Please visit our website at: www.bpsbioscience.com