

Data Sheet

TGF β R2

Human, recombinant, N-terminal GST-tag
Catalog #: 40707
Lot#: 140205-G1 Conc.: 0.11 mg/ml

Formulated in: 40 mM Tris pH 8, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 3 mM DTT, and 20% glycerol

Stability: >6 months at -80°C . Avoid freeze/thaw cycles.

References:

1. Lee J, *et al.*, *PLoS ONE*. 2013 **8(2)**:e57074.
2. Bhowmick, N.A, *et al.*, *Science*. 2004; **303(5659)**:848-851

Description:

Human transforming growth factor beta receptor 2 (TGF β R2), also known as TGFBR2, TGFR-2, and TGF-Beta Receptor Type 2 GenBank Accession No. NM_003242 (var2), a.a. 190-567 (end), with N-terminal GST-tag, MW=69.9 kDa, expressed in an Sf9 cell expression system.

Application:

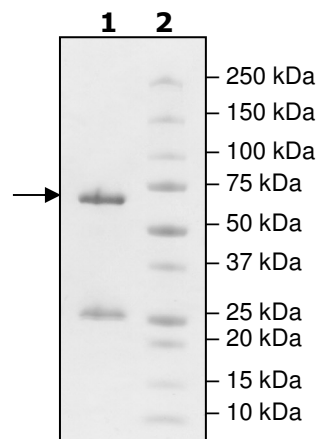
Useful for studying protein binding, receptor activity, and screening small molecules.

Quality Assurance

4-20% SDS-PAGE

Lane 1:
2.2 μg TGF β R2
Lane 2:
Protein Marker

MW: 69.9 kDa
Purity: $\geq 66\%$



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



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

Protein Sequence:

MSPILGYWKIKGLVQPTRLLEYLEEKYEEHLYERDEGDKWRNKKFELGLEFPN
LPYYIDGDVKLTQSMAIRYIADKHNMLGGCPKERAEISMLEGAVLDIRYGVSRIA
YSKDFETLKVDFLSKLPEMLKMFKDRLCHKTYLNGDHVTHPDFMLYDALDVVLY
MDPMCLDAFPKLVCFKKRIEAIPQIDKYLKSSKYIAWPLQGWAQATFGGGDHLEV
LFQGPLGSRVNRQQKLSSTWETGKTRKLMFSEHCAIILED RSDISSTCANNI
NHNTELLPIELDTLVGKGRFAEVYKAKLKQNTSEQFETVAVKIFPYEYASWK
TEKDIFSDINLKHENILQFLTAERKTEL GKQYWLITAFHAKGNLQEYLTRHVIS
WEDLRKLGSSLARGIAHLHSDHTPCGRPKMPIVHRDLKSSNILVKNDLTCCLC
DFGLSLRLDPTLSVDDLANSQVGTARYMAPEVLES RMNLENVESFKQTDVY
SMALVLWEMTSRCNAVGEVKDYEPFSGSKVREHPCVESMKDNVLRDRGRPE
IPSFVLNHQGIQMCETLTECWDHDPEARLTAQCVAERFSELEHLDRLSGRS
CSEKIPEDGSLNTTK

Black: TGF β R2 (a.a. 190-567)

Red: Flag Fusion

Blue: Linker including PreScission Protease Cleavage Site

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