

Data Sheet

PDGFR α (D842I), GST-Tag

Human, Recombinant, N-terminal GST-Tag
Catalog #: 100202
Lot #: 181112-G1 **Conc.:** 0.33 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0,
110 mM NaCl, 2.2 mM KCl, 0.04% Tween-
20, 3mM DTT and 20% glycerol.

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

Description: Recombinant human PDGFR α , also known as PDGFR2, RHEPDGFRA, Platelet-derived growth factor receptor alpha, and CD140a, with D842I mutation, Genbank accession number NM_006206, a.a. 550-1089 (end), with an N-terminal GST-tag, expressed in an Sf9 insect cells via a baculovirus expression system. MW = 88 kDa.

References:

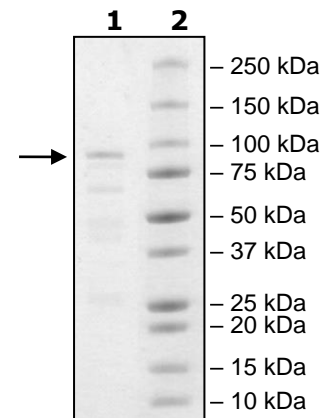
1. Calabuig-Fariñas, S. *et al. Virchows Archiv.* 2009 Mar; **454(3)**:327-331
2. Hojjat-Farsangi, M. *Int. J. Mol. Sci.* 2014 Aug; **15(8)**:13768-13801

Quality Assurance

4-20% SDS-PAGE Coomassie staining

Lane 1:
2 μg PDGFR α
Lane 2:
Protein Marker

MW: 88 kDa
Purity: $\geq 60\%$



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

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

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

Please visit our website at: www.bpsbioscience.com