

## Data Sheet

### PDGFR $\alpha$ (D842Y), GST-Tag

Human, Recombinant, N-terminal GST-tag  
**Catalog #:** 100201  
**Lot #:** 181030-G      **Conc.:** 0.15 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  $-80^{\circ}\text{C}$ . Avoid  
 freeze/thaw cycles.

**References:**

1. Floege, J. *et al. J Am Soc Nephrol.* 1998  
 Feb; **9(2)**:211-23.
2. Kumagai, S. *et al. Tohoku J Exp Med.*  
 2001 Sep; **195(1)**:21-33.

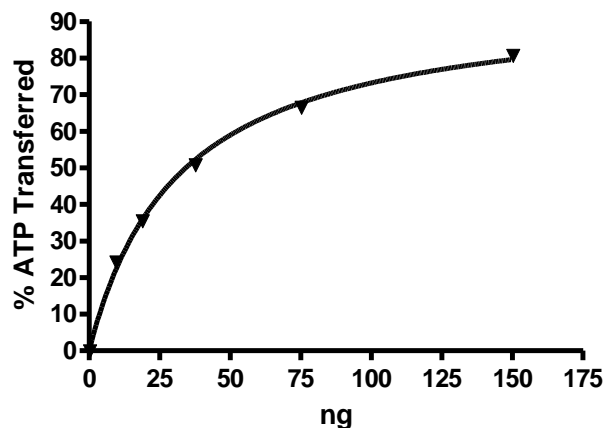
**Description:** Recombinant human  
 PDGFR $\alpha$ , also known as CD140A, and  
 PDGFR2, with D842Y mutation,  
 Genbank accession number  
 NM\_00626, with an N-terminal GST-  
 tag, expressed in Sf9 insect cells via a  
 baculovirus expression system. MW=88  
 kDa.

**Specific Activity:** 153 pmole/min/ $\mu\text{g}$   
 Assay Conditions: Assay was done in  
 Kinase buffer #79334 using Poly-  
 (Glu4:Tyr) substrate (0.2 mg/ml) and 10  
 $\mu\text{M}$  ATP. Reaction was done at  $30^{\circ}\text{C}$  for  
 45 min. Amount of ATP transferred was  
 calculated using 50  $\mu\text{l}$  Kinase-Glo  
 reagent (Promega).

**Applications:** Useful for the study of  
 enzyme kinetics, screening inhibitors,  
 and selectivity profiling.

## Quality Assurance

### Specific Activity

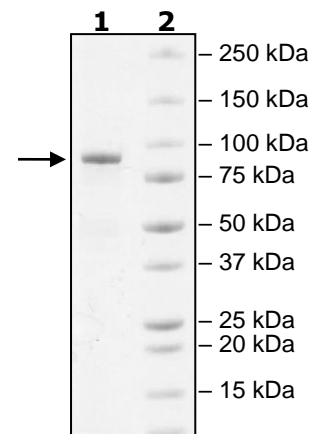


### 4-20% SDS-PAGE Coomassie staining

**Lane 1:**  
 2 $\mu\text{g}$  PDGFR $\alpha$   
**Lane 2:**  
 Protein Marker

**MW:** 88 kDa

**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 **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)