

## Data Sheet

### p120GAP, His-Tag

Human, Recombinant, N-terminal His-Tag

**Catalog #:** 100518

**Lot#:** 200102-1

**Conc.:** 0.21 mg/ml

**Formulated in:** 40 mM Tris-HCl pH 8.0, 110 mM NaCl, 2.2 mM KCl, 20% glycerol.

**Stability:** At least 6 months at  $-80^{\circ}\text{C}$ .  
 Avoid freeze/thaw cycles.

**References:**

1. Scheffzek, K., *et al.*, *Nature*. 1996; **384**: 591-596
2. Torit, M. *et al.*, *JBC*. 1992; **267**: 8293-8298

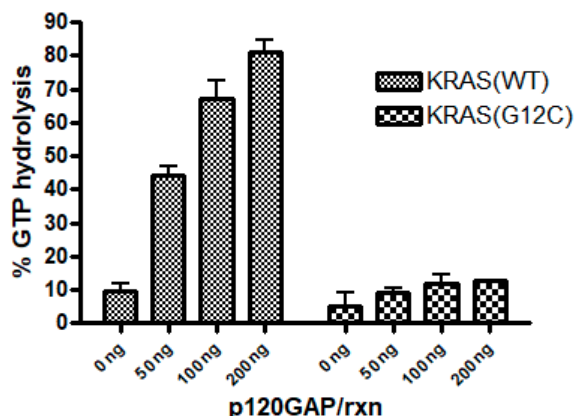
**Description:** Human Ras GTPase-activating protein 1, also known as p120GAP, RASA1, GAP and RASA, GenBank Accession No. NM\_002890, a.a. 714-1047(end) with an N-terminal His-Tag, expressed in an *E. coli* cell expression system. MW=39 kDa.

**Assay Conditions:** GTPase activity of KRAS(WT, 50ng) and KRAS(G12C, 100ng) were tested in the presence of various concentration of p120GAP (0 ng - 200 ng) and 2  $\mu\text{M}$  GTP by using GTPase-Glo reagents (Promega). Reactions were done at room temperature for 80 min.

**Applications:** Useful for the study of RAS GTPase activation.

### Quality Assurance

**GAP assisted GTPase activity**

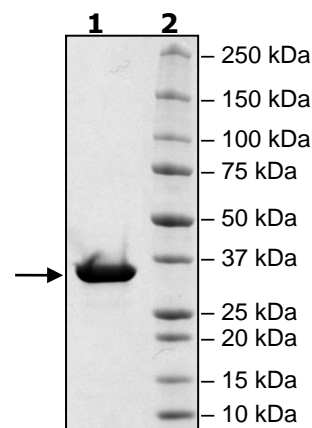


**4-20% SDS-PAGE**  
**Coomassie staining**

**Lane 1:**  
 2 $\mu\text{g}$  p120GAP  
**Lane 2:**  
 Protein Marker

**MW:** 39 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)