

Data Sheet

NRAS (Q61K), His-Tag

Human, Recombinant, N-terminal His-tag

Catalog #: 100800

Lot #: 200813

Conc: 2.35 mg/ml

Description: Human NRAS also known as GTPase NRas and NS6, GenBank Accession No. NM_002524, a.a. 2-186, with Q61K mutation and N-terminal His-tag, expressed in a Baculovirus infected Sf9 cell expression system. MW=22 kDa.

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween20, 20% glycerol, 3 mM DTT.

Stability: At least 6 months at -70°C. Avoid freeze/thaw cycles.

References:

1. Roock, W. D. *et al.*, *The Lancet Oncology*. 2020 Aug; **11(8)**: 753-762.
2. Kumari, S. *et al.*, *Nat. Chem. Bio.* 2007 Feb; **3**: 218-221.

Quality Assurance

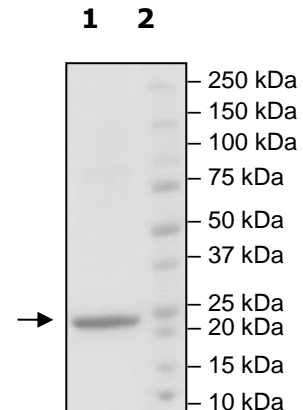
6-20% SDS-PAGE Coomassie staining

Lane 1:
NRAS (Q61K)

Lane 2:
Protein Marker

Purity: >90%

MW: 22 kDa



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