

Data Sheet

DRAK2

Human, Recombinant

Catalog #: 71225

Lot #: 160804A1

Conc: 0.5 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 20% glycerol, 1 mM DTT, 1 mM EDTA
Stability: At least 6 months at -80°C . Avoid freeze/thaw cycles. Storing diluted enzyme is not recommended. If necessary, use carrier protein (BSA 0.1-0.5%).

References:

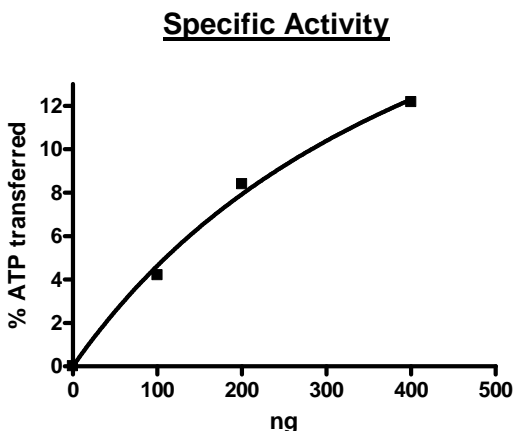
1. McGargill, M. *et al.*, *Immunity* **21** (6): 781-791 (2004).
2. Mao, J. *et al.*, *J. Biol. Chem.* **281**: 12587-12595 (2006).

Description: Human DRAK2, also known as DAP kinase-related apoptosis-inducing protein kinase 2 or death-associated protein kinase-related 2 (GenBank Accession No. NM_004226.3), a.a. 2-372 (end), MW = 43 kDa, expressed in a baculovirus-infected Sf9 cell expression system.

Assay Description: Assay was done in Kinase buffer containing 1 mM DTT using MRCL3 peptide substrate (0.2 mg/ml) and 20 μM ATP. Reaction was done at 30°C for 50 min. Amount of ATP transferred was calculated using Kinase-Glo reagent (Promega)

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

Quality Assurance



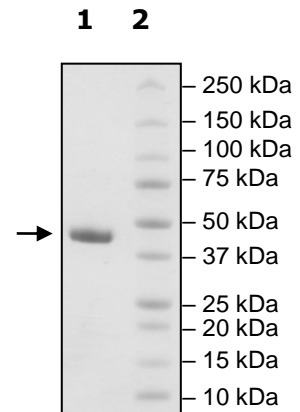
4-20% SDS-PAGE
Coomassie staining

Lane 1:
DRAK2

Lane 2:
Protein Marker

Purity: $\geq 90\%$

MW: 43 kDa



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

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

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

Please visit our website at: www.bpsbioscience.com