

Data Sheet

MRCK α , GST-tag

Human, Recombinant, with N-terminal GST-tag

Catalog #: 40129

Lot #: 170620 **Conc.:** 0.1 mg/ml

Formulated in: 50 mM Tris-HCl, pH 7.5, 150 mM NaCl, 0.25 mM DTT, 0.1 mM EGTA, 0.1 mM EDTA, 0.1 mM PMSF, 25% glycerol.

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:

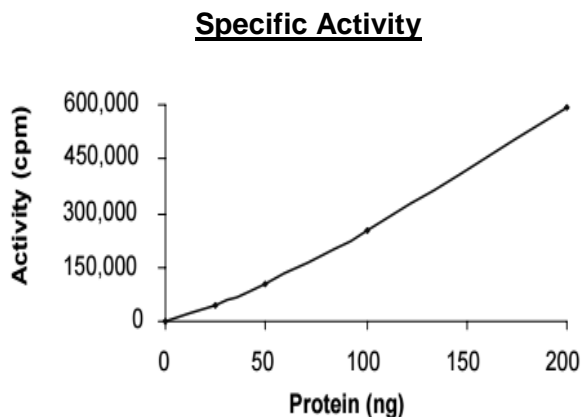
1. Ivan, T. et al: Gene; 2003; 304:107-115.
2. Ivan, T. et al. J. Biol. Chem., 2001: 276 (24): 21209-21216.

Description: Recombinant human Myotonic Dystrophy Kinase-Related Cdc42-Binding Kinase alpha (MRCK α), a.a. 1-473, with N-terminal GST-tag, expressed in Sf9 insect cells via a baculovirus expression system.

Specific Activity: 101 pmol/min/ μg .

Application: Useful for studying enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Assurance

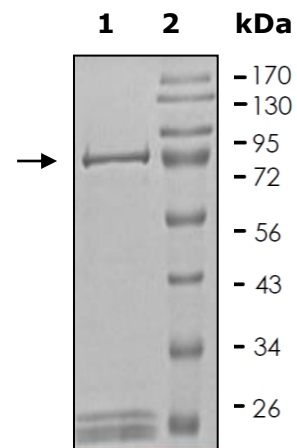


10% SDS-PAGE Coomassie staining

Lane 1:
MRCK alpha

Lane 2:
Protein Marker

MW: 73 kDa
Purity: $\geq 75\%$



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