

Data Sheet

DRAK2, GST-Tag

Human, Recombinant, N-terminal GST-tag
Catalog #: 71224
Lot #: 190530-4 **Conc:** 0.62 mg/ml

Formulated in: 50 mM HEPES, pH 7.5, 0.001% Brij-35, 1 mM DTT, 10 mM MgCl₂, 150 mM KCl, 5% Glycerol, 0.1 mM EDTA.

Stability: At least 6 months at -80°C. Avoid freeze/thaw cycles.

References:

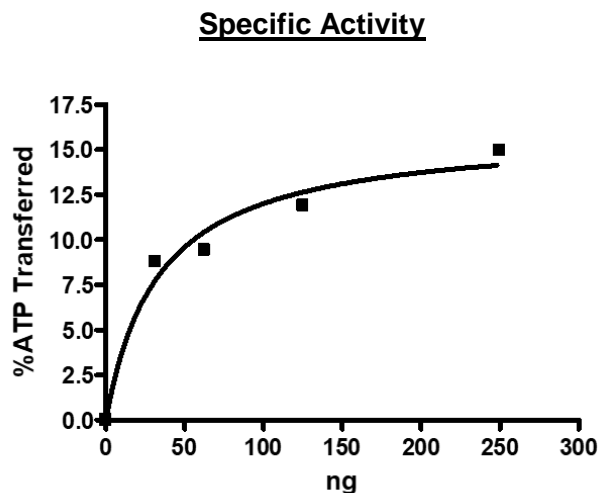
1. McGargill, M. et al, Immunity. **21 (6):** 781-791 (2004).
2. Mao, J. et al, Journal of Biological Chemistry. **281:** 12587-12595 (2006).

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), with N-terminal GST-tag, expressed in a baculovirus-infected Sf9 cell expression system. MW = 69 kDa.

Specific Activity: 15 pmole/min/μg
Assay Conditions: Assay was done in Kinase buffer containing 1 mM DTT using MRCL3 peptide substrate (0.2 mg/ml) 20 μM ATP. Reaction was done at 30°C for 45 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:
2 μg DRAK2
Lane 2:
Protein Marker
MW: 69 kDa
Purity: 75%

