

PRODUCT DATA SHEET



PAK7

CATALOG NO.: SE-390

LOT NO.: P8615

DESCRIPTION: PAK7, MW=104 kDa, full-length human recombinant containing an N-terminal GST tag expressed in insect cells. GenBank ID BC024179.

PURITY: >75% by SDS-PAGE.

SPECIFIC ACTIVITY: 55 nmol/min/mg. Kinase activity was measured as the molar amount of phosphate incorporated into a peptide substrate (CKRPRAASFAE) at 30°C using a concentration of 50 μ M [³²P]ATP in 25 mM MOPS, pH 7.2, 12.5 mM β -glycerophosphate, 5 mM EGTA, 2 mM EDTA, 25 mM MgCl₂, 0.25 mM DTT (added right before using).

Deleted: 4 mM MOPS, pH 7.2, 2.5 mM β -glycerophosphate, 1 mM EGTA, 0.4 mM EDTA, 4 mM MgCl₂, 0.05 mM DTT.

APPLICATIONS: Study enzyme regulation and kinetics, phosphorylate target substrates, screen for inhibitors.

SUPPLIED AS: 0.1 mg/ml 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 and 25% glycerol.

STORAGE: -80°C. After initial defrost, aliquot product into individual tubes and refreeze the remaining, unused enzyme quickly by snap-freezing in a dry/ice ethanol bath or liquid nitrogen, if possible. Avoid repeated freeze/defrost cycles.

NOTE: When stored under the above conditions, this enzyme is stable at the concentration supplied, in its current storage buffer. Procedures such as dilution of the enzyme followed by refreezing, could lead to loss of activity.

FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.