

Data Sheet

PKC α , GST-Tag

Human, Recombinant, N-terminal GST tag
Catalog #: 40157
Lot#: 170329-3 **Conc.:** 0.52 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween20, 3 mM DTT, 20% glycerol, 1.6 mM Glutathione

Stability: At least 6 months at -80°C . Avoid freeze/thaw cycles. Protein may be diluted to $\geq 100 \mu\text{g/ml}$ in PBS + glycerol and stored at -80°C .

References:

1. Braz, J. Et al., Nature Med. 10:248-254 (2004).
2. Coussens, L. et al., Science.233:859-866 (1986).

Description:

Recombinant human PKC α , full length, (GenBank Accession No. NM_002737) with N-terminal GST-tag, expressed in Sf9 insect cells via a baculovirus expression system. MW= 102 kDa.

Specific Activity: $\geq 250 \text{ pmol/min}/\mu\text{g}$

One unit will transfer 1pmol of phosphate from ATP to Histone H1 in 1 minute at pH 7.4 at 30 degrees C.

250units/ug protein

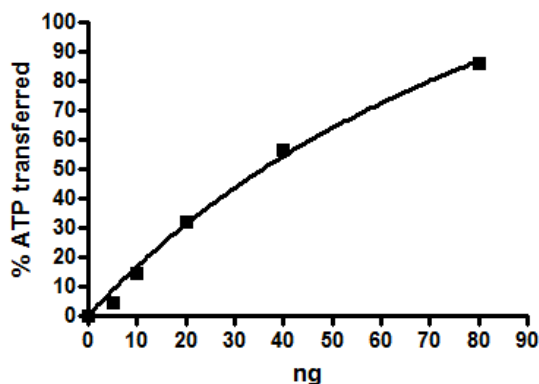
Assay Conditions: Assay was done in Kinase buffer containing 0.1 mM CaCl₂ and lipid activator (1 mg/ml Phosphatidyl Serine and 0.02 mg/ml Diacylglycerol in 0.3% TritonX-100) using Histone H1 (0.1 mg/ml) and 20 μM ATP. Reaction was done at 30°C for 40 min. Amount of ATP transferred was calculated using Kinase-Glo reagent (Promega)

Application:

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

Quality Assurance

Specific Activity

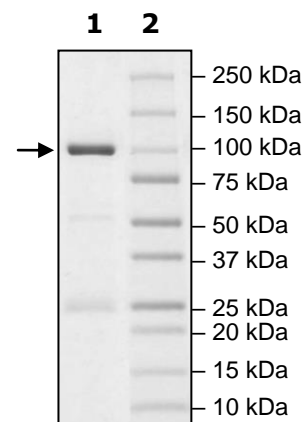


4-20% SDS-PAGE Coomassie staining

Lane 1: PKC α
Lane 2: Protein Marker

Purity: $\geq 80\%$

MW: 102 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