

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.

References:

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

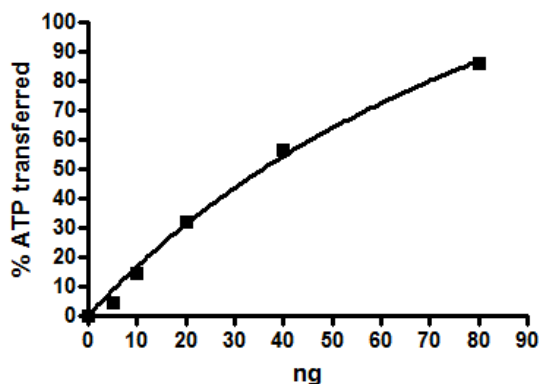
Description: Human PKC α , GenBank
 Accession No. NM_002737, full length,
 with N-terminal GST-tag, expressed in
 Sf9 insect cells via a baculovirus
 expression system. MW= 102 kDa.

Specific Activity: ≥ 320 pmol/min/ μg
 One unit will transfer 1pmol of phosphate
 from ATP to Histone H1 in 1 minute at
 pH 7.4 at 30°C . 250 units/ μg 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).

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

Quality Assurance

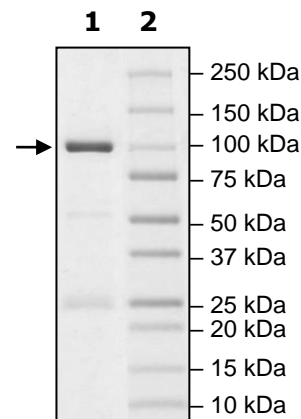
Specific Activity



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 **1.858.202.1401** Fax **1.858.481.8694**
 Or you can Email us at: info@bpsbioscience.com
 Please visit our website at: www.bpsbioscience.com