

Data Sheet

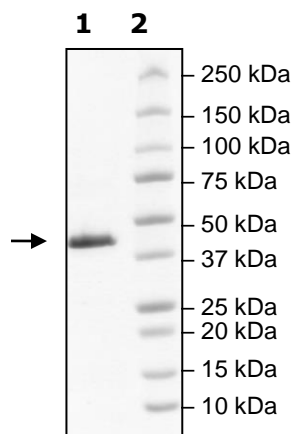
PKA α , His-Tag

Human, Recombinant, N-terminal His-Tag

Catalog #: 79044**Lot #:** 170601**Conc.:** 0.98mg/ml**Description:** Human PKA α , GenBank Accession No. NM_002730, 2-351 (end), with N-terminal His-Tag, MW = 42 kDa, expressed in an E.coli expression system.**Formulated in:** 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween20, 3mM DTT, 20% glycerol, 64mM imidazole**Stability:** At least 6 months -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. Tasken, K. et al, 19p13.1. Genomics 36: 535-538, 1996.
2. Skalhegg, BS. et al, Molec. Endocr. 16: 630-639, 2002.

Quality Assurance

**4-20% SDS-PAGE
Coomassie staining****Lane 1:**
PKACA, WT**Lane 2:**
Protein Marker**Purity:** $\geq 90\%$ **MW:** 42 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.comPlease visit our website at: www.bpsbioscience.com



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Sequence

MASRGSHHHHHHGS MGNAAAAKKGSEQESVKEFLAKAKEDFLKKWESPAQNTA
HLDQFERIKTLGTGSFGRVMLVKHKETGNHYAMKILDKQKVVKLKQIEHTLNE
KRILQAVNFPFLVKLEFSFKDNSNLYMVMEYVPGGEMFSHLRRIGRFSEPHAR
FYAAQIVLTFEYLSLDLIYRDLKPENLLIDQQGYIQVTDFGFAKRVKGRWT
LCGTPEYLAPEIILSKGYNKAVDWWALGVLIYEMAAGYPPFFADQPIQIYEKI
VSGKVRFP SHFSSDLKDLLRNLLQVDLTKRFGNLKNGVNDIKNHKWFATTDWI
AIYQRKVEAPFIPKFKGPGDTSNFDYEEEEIRVSINEKCGKEFSEFLEA

Red: His-tag and a short linker

Black: PKACA (aa 2-351(end))

S

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