

# PKC-theta

aa1-706

**Product No.:** 0113-0000-1

**Lot:** 008

**Description:** Human PKC-theta

Amino acids M<sub>1</sub>-S<sub>706</sub> (as in GenBank entry NM\_006257)\*, N-terminally fused to GST-FactorXa cleavage site, C-terminally fused to His<sub>6</sub>

\*Sequence may contain documented polymorphisms  
Detailed sequence on request

**Theoretical MW<sub>Fusion Protein</sub>:** 109,677 Da

**Product identity:** PKC-theta, Lot 008, was confirmed as human PKC-theta by mass spectroscopy LC-ESI-MS/MS (Protagen AG, Germany)

**Expression:** Baculovirus infected Sf9 cells

**Purification:** One-step affinity purification using GSH-agarose

**Storage buffer:** 50 mM Tris-HCl, pH 8.0; 100 mM NaCl, 5 mM DTT, 4 mM reduced glutathione, 20% glycerol

**Storage temperature:** -80°C  
Avoid repeated freeze-thaw cycles!

**Protein concentration:** 0.273 µg/µl  
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

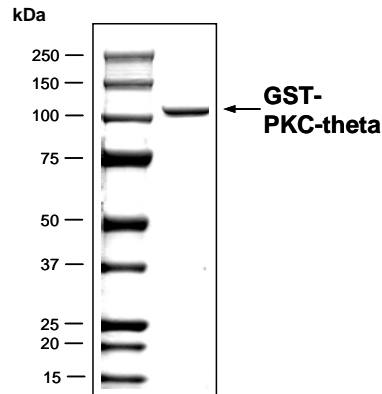
**Method for determination of K<sub>m</sub> value & Specific activity:**

- Assay conditions:
  - 60 mM HEPES-NaOH, pH 7.5
  - 3 mM MgCl<sub>2</sub>
  - 3 mM MnCl<sub>2</sub>
  - 1 mM CaCl<sub>2</sub>
  - 4 mM EDTA
  - 3 µM Na-orthovanadate
  - 5 µg / ml Phosphatidylserine
  - 1 µg / ml 1.2 Dioleoyl-glycerol
  - 1.2 mM DTT
  - 50 µg / ml PEG<sub>20,000</sub>
  - ATP (variable)
  - Substrate: Histone H2B, 30 µg / ml
  - Recombinant PKC-theta: 4.0 µg / ml

- Filter binding assay  
MSFC membrane (Millipore)

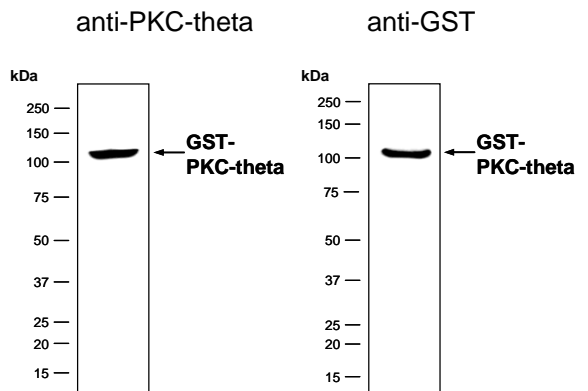
**Specific activity:** 166 pmol/µg×min

**Coomassie stain:**



2.0 µg GST-PKC-theta

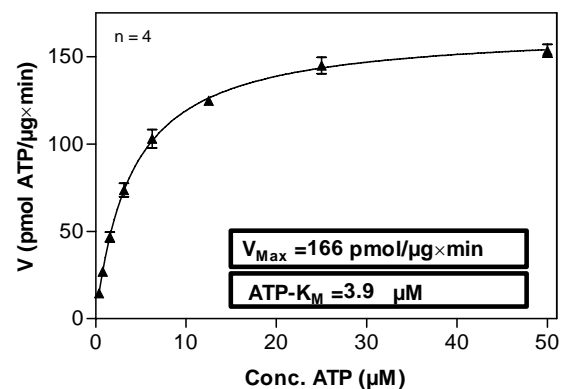
**Western blot analysis:**



500 ng GST-PKC-theta

500 ng GST-PKC-theta

**Determination of K<sub>m</sub> value for ATP:**



This product is for in vitro research use only, not for use in humans or animals. ProQinase disclaims any warranty explicitly or implied that the use of the product or parts of the product is free from third party intellectual property claims unless this is explicitly stated.