

Aurora-A

aa1-403

Product No.: 0166-0000-1

Lot: 004

Description: Human Aurora-A
Amino acids M₁-S₄₀₃ (as in GenBank entry NM_003600)*, N-terminally fused to GST-HIS₆-Thrombin cleavage site

*Sequence may contain documented polymorphisms
Detailed sequence on request

Product identity: Aurora-A, Lot 004, was confirmed as Aurora-A by specific Western Blotting using anti Aurora-A antibody

Theoretical MW_{Fusion Protein}: 76,334 Da

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, 15 mM reduced glutathione, 20% glycerol

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

Protein concentration: 0.245 µg/µl
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

Method for determination of K_m value & Specific activity:

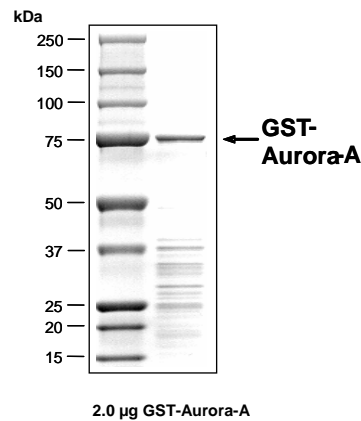
• Assay conditions:

60 mM HEPES-NaOH, pH 7.5
3 mM MgCl₂
3 mM MnCl₂
3 µM Na-orthovanadate
1.2 mM DTT
2.5 µg / 50 µl PEG_{20,000}
ATP (variable)
Substrate: tetra(LRRLSLG), 5 µg / 50 µl
Recombinant Aurora-A: 200 ng / 50 µl

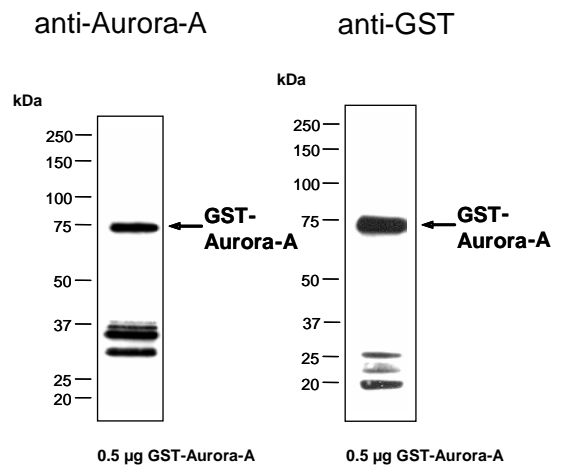
• Filter binding assay
MAPH membrane (Millipore)

Specific activity: 129 pmol/µg×min

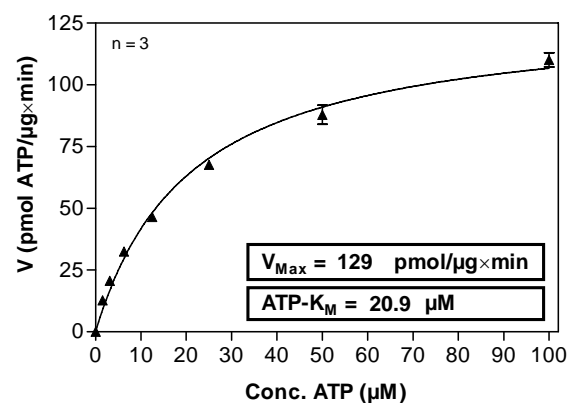
Coomassie stain:



Western blot analysis:



Determination of K_m 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.