

Certificate of Analysis



ATF2 (activating transcription factor 2)

Recombinant Protein Kinase Substrate

Synonyms: CRE-BP1, CREB2, TREB7

Product No.: 0594-0000-2

Lot: 010

Description: Human ATF2, N-terminal fragment, amino acids M₁₉-D₁₁₀ (as in GenBank entry NM_001880)*, N-terminal GST-HIS₆ fusion protein with a 3C cleavage site, expressed in E.coli

*Sequence may contain documented polymorphisms
Detailed aa-sequence on request

Theoretical MW: 38,620 Da

Expression: E.coli

Purification: Affinity purification using GSH-agarose

ATPase activity: In an ADP-Glo™ assay (Promega) with 10 μM ATP or 30 μM ATP, the ATP → ADP conversion within 30 min is below 1% at a concentration of 100 μg/ml ATF2*

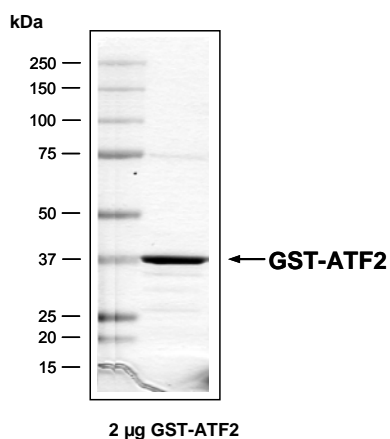
*detailed ATPase assay conditions on request

Storage buffer: 50 mM HEPES pH 7.5, 100 mM NaCl, 1 mM DTT, 10% glycerol, 15 mM GSH

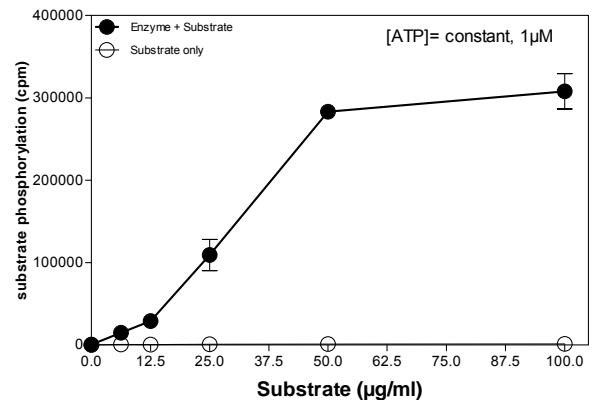
Storage temperature: -80°C
Avoid repeated freeze-thaw cycles!

Protein concentration: 0.648 μg/μl
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

Coomassie stain:



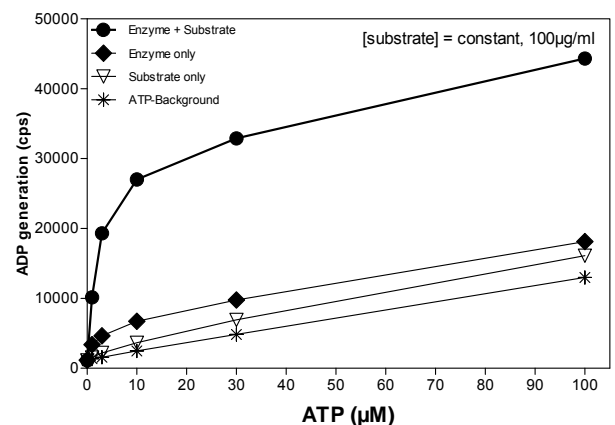
Phosphorylation of ATF2 by the kinase p38-beta (Radiometric filter binding assay):



Assay mixture:

70 mM HEPES-NaOH, pH 7.5
3 mM MgCl₂
3 mM MnCl₂
3 μM Na-orthovanadate
1.2 mM DTT
50 μg/ml PEG_{20,000}
ATP: 1 μM
Substrate (ATF2): variable concentration
p38-beta: 0.5 μg/ml
MSFC membrane (Millipore)

Phosphorylation of ATF2 by the kinase p38-beta (ADP-Glo™ assay / Promega):



Assay mixture:

70 mM HEPES-NaOH, pH 7.5
3 mM MgCl₂
3 mM MnCl₂
3 μM Na-orthovanadate
1.2 mM DTT
50 μg/ml PEG_{20,000}
ATP: variable concentration
1 % (v/v) DMSO
Substrate (ATF2): 100 μg/ml
p38-beta: 1 μg/ml

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.