

## ProQinase™ ATF2

activating transcription factor 2

### Recombinant Protein Kinase Substrate

**HGNC Symbol:** ATF2

**Synonyms:** CRE-BP1, CREB2, HB16, TREB7

**Product No.:** 0594-0000-2

**Lot:** 014

**Description:** Human ATF2, N-terminal fragment, amino acids M<sub>19</sub>-D<sub>110</sub> (as in [NCBI/Protein](#) entry NP\_001871.2), N-terminal GST-HIS<sub>6</sub> fusion protein with a 3C cleavage site, expressed in E.coli

**Theoretical MW<sub>Fusion Protein</sub>:** 38,620 Da

**Expression host:** E.coli

**Purification:** GST-Affinity Chromatography

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

Detailed ATPase assay conditions on request

**Storage buffer:** 50 mM HEPES pH 7.5, 100 mM NaCl, 5 mM DTT, 15 mM reduced glutathione, 10 % glycerol

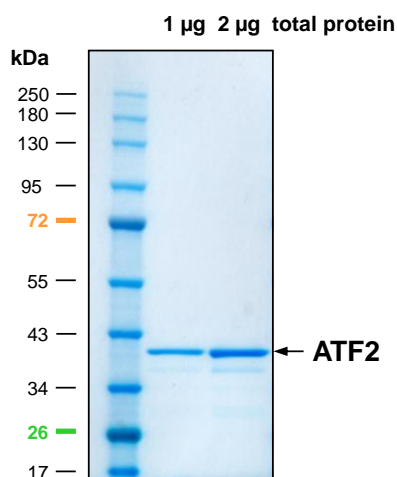
**Storage temperature:** -80°C

For complete recovery, mix well and spin before use. Product must not be stored in diluted solutions, aliquots below 10 μl are not advisable. Avoid repeated freeze-thaw cycles!

**Protein concentration:** 2.25 μg/μl  
(Bradford method using BSA [Sigma, cat# A-7638, Lot 79H7641] as standard protein)

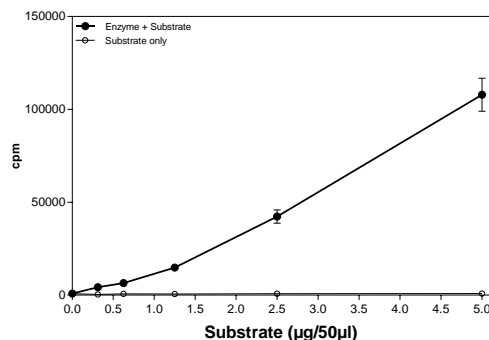
**ATF2 Lot 014:**

**Coomassie stain**



### Phosphorylation of ATF2 Lot014 by p38beta

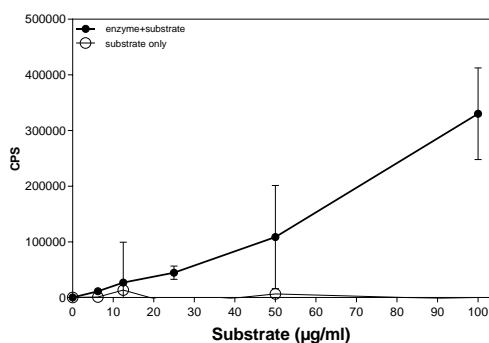
#### Radiometric filter binding assay



#### Assay conditions:

70 mM HEPES-NaOH, pH 7.5  
3 mM MgCl<sub>2</sub>  
3 mM MnCl<sub>2</sub>  
3 μM Na-orthovanadate  
1.2 mM DTT  
50 μg/ml PEG<sub>20,000</sub>  
ATP: 1 μM  
Substrate: variable concentration  
Kinase: 1 μg/ml  
MSFC filter plate

#### ADP-Glo™ assay (Promega)



#### Assay conditions:

70 mM HEPES-NaOH, pH 7.5  
3 mM MgCl<sub>2</sub>  
3 mM MnCl<sub>2</sub>  
3 μM Na-orthovanadate  
1.2 mM DTT  
50 μg/ml PEG<sub>20,000</sub>  
ATP: variable concentration  
1 % (v/v) DMSO  
Substrate: 100 μg/ml  
Kinase: 1 μg/ml

# ATF2

Product No.: 0594-0000-2

ATF2 Recombinant Fusion Protein Amino Acid Sequence							
1	MSPILGYWKI	KGLVQPTRL	LEYLEEKYEE	HLYERDEGDK	WRNKKFELGL	EFPNLPYYID	60
61	GDVKLTQSMA	IIRYIADKHN	MLGGCPKERA	EISMLEGAVL	DIRYGVSRIA	YSKDFETLKV	120
121	DFLSKLP EML	KMFEDRLCHK	TYLNGDHVTH	PDFMLYDALD	VVLYMDPMCL	DAFPKLVCFK	180
181	KRIEAI PQID	KYLKSSKYIA	WPLQG WQATF	GGGDHPPKSD	PMGHHHHH G	RDSLEVL FQG	240
241	PLAMSDDKPF	LCTAPGCGQR	FTNEDHLAVH	KHKHEMTLKF	GPARNDSVIV	ADQTPTPTRF	300
301	LKNCEEVGLF	NELASPFENE	FKKASEDDIK	KMPLD			360

1-218: GST   **Red**: HIS6-tag   **Green**: 3C cleavage site   **blue**: ATF2 fragment

This product was manufactured at Reaction Biology in Freiburg, Germany, and is for in vitro research use only, not for use in human or animals.  
 © European Union, 2020. Material may not be reproduced or distributed without written permission from Reaction Biology Europe GmbH