

## Data Sheet

### ATF2

Human, recombinant, N-terminal GST tag

**Catalog #:** 40520

Lot #: 120716D      Conc.: 0.88 mg/ml

**Formulated in:** 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 16 mM Glutathione, 3 mM DTT, and 20% Glycerol.

**Stability:** >6 months at -80°C

#### References:

1. Huang, Y., *et al.*, *J Mol Signal.* 2008. **12**:3:3.
2. Maekawa, T., *et al.*, *Oncogene.* 2008. **27**(8):1045-54.
3. Clerk, A. And Sugden, P.H. *Biochem. J.* 1997. **325**: 801-810.

#### Description:

Human ATF2 (GenBank Accession No. NM\_001880), (amino-acids 19 - 96) with N-terminal GST tag, MW= 34.8 kDa, expressed in *E. coli* expression system.

#### Application:

A suitable substrate for JNK1 and JNK2.

ATF2 has been shown in the literature to appear as multiple bands due to differential phosphorylation.<sup>3</sup>

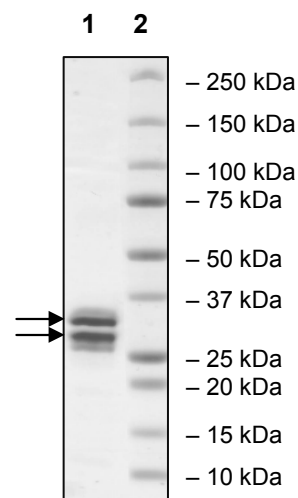
## Quality Assurance

### 4-20% Gradient SDS-PAGE Coomassie staining

**Lane 1:**  
2 µg ATF2

**Lane 2:**  
Protein Marker  
Biorad (#161-0373)

**MW:** 34.8 kDa  
**Purity:** ≥85%



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](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)