

Product Information

logo

ATF7 Transcription Factor Monoclonal Antibody

Cat No: G01001 - 100 µL

General Data

Shipping:	wet ice
Stability:	Store at -20degrees; shelf life 3 year maximum after production Avoid repeat freezing and thawing cycles. Does not contain any preservative
Formulation:	Liquid Ascite, does not contain any preservative therefore avoid repeat freezing and thawing cycles
Host:	Mouse
Antigen:	recombinant human ATF7 protein
Clone:	2F10
Isotype:	IgG1, κ
Application(s):	ELISA Western Blot Immunofluorescence Immunoprecipitation gel supershift recommended dilutions: 1/500-1/5000
Specificity:	human ATF7 protein, epitope localized between aa 257 and 398

Product Overview

The activating transcription factors (ATFs) belong to the AP-1 family of transcription factors. ATF consists of seven members, ATF1 to 7. ATF7 (originally called ATFa) is ubiquitously expressed in various tissues and plays important roles in G1 and S phases. ATF7 might be important in controlling memory in cells of the innate immune system.

FP/18/20

For research laboratory use only - Not for human diagnostic use.

Buyers agree to purchase the material subject to Purchasing Terms that can be found on our website. Seek appropriate training to safely handle this product under normal conditions. Use the recommended personal protective equipment to prevent chemical exposures.

Bertin Bioreagent does not make any other warranty or representation whatsoever whether expressed or implied, with respect to these products. In no event will Bertin Bioreagent be liable for incidental, consequential or punitive damages.

Contact Bertin Bioreagent

Parc d'Activités du Pas du Lac
10 bis avenue Ampère
78180 Montigny le Bretonneux - France
Tel.: +33 (0)139 306 036
<https://www.bertin-bioreagent.com/pa206/contact-us>
<https://www.bertin-bioreagent.com>

Product Information

logo

ATF7 Transcription Factor Monoclonal Antibody

Cat No: G01001 - 100 µL

Scientific Literature

Hamard PJ, Dalbies-Tran R, Hauss C, Davidson I, Kedinger C, Chatton B. A functional interaction between ATF7 and TAF12 that is modulated by TAF4. *Oncogene*. 2005 May 12;24(21):3472-83.

FP/18/20

For research laboratory use only - Not for human diagnostic use.

Buyers agree to purchase the material subject to Purchasing Terms that can be found on our website. Seek appropriate training to safely handle this product under normal conditions. Use the recommended personal protective equipment to prevent chemical exposures.

Bertin Bioreagent does not make any other warranty or representation whatsoever whether expressed or implied, with respect to these products. In no event will Bertin Bioreagent be liable for incidental, consequential or punitive damages.

Contact Bertin Bioreagent

Parc d'Activités du Pas du Lac
10 bis avenue Ampère
78180 Montigny le Bretonneux - France
Tel.: +33 (0)139 306 036
<https://www.bertin-bioreagent.com/pa206/contact-us>
<https://www.bertin-bioreagent.com>