



TAFII 68Å Monoclonal Antibody

Cat No: G01089 - 100 ÅµL

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:	Synthetic peptide coupled to ovalbumin
Clone:	2B11
Isotype:	IgG1, κ



Application(s):	ELISA Western Blot (endogenous and recombinant proteins) Immunocytofluorescence Immunoprecipitation Recommended dilutions: 1/500-1/5000
------------------------	---

Specificity: Human, other species have not been tested

Scientific Literature

Bertolotti, A., Lutz, Y., Chambon, P. and Tora, L. (1996) EMBO J. 15, 5022-5031.

Bertolotti A., Melot T., Acker J., Vigneron M., Delattre O. and Tora L. (1998) Mol. Cell. Biol. 18, 1489-1497.

FP/03/20

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>