



TAF11^Δ Monoclonal Antibody

Cat No: G01077 - 100 μ L

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 TAF11 protein
Clone:	2B4
Isotype:	IgG1, κ



Application(s):	ELISA
	Western Blot
	Immunoprecipitation
	Immunofluorescence
	Recommended dilutions: 1/500-1/5000

Specificity: human and mouse TAF11

Scientific Literature

Mengus G., May M., Jacq X., Staub A., Tora L., Chambon P. and Davidson I. (1995). Cloning and characterization of hTAFII18, hTAFII20 and hTAFII28: three subunits of the human transcription factor TFIID. EMBO J., 14, 1520-1531

May M., Mengus G., Lavigne A.

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>