



TAF4 (formerly TAFII135) Monoclonal Antibody

Cat No: G01083 - 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 TAF4 protein
Clone:	1B12
Isotype:	IgG1, κ



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

Specificity: human and mouse TAF4

Scientific Literature

Mengus G, May M, Carre L, Chambon P, Davidson I. Human TAF II 135 potentiates transcriptional activation by the AF-2s of the retinoic acid, vitamin D3, and thyroid hormone receptors in mammalian cells. *Genes Dev.* 1997, 11, 1381-95.

Mazzarelli JM, Mengus G

FP/10/20

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>