



## TIF2 Monoclonal Antibody

Cat No: G01104 - 100 µL

### General Data

<b>Shipping:</b>	wet ice
<b>Stability:</b>	Store at -20degrees; shelf life 3 year maximum after production Avoid repeat freezing and thawing cycles. Does not contain any preservative
<b>Formulation:</b>	Liquid Ascite, does not contain any preservative therefore avoid repeat freezing and thawing cycles
<b>Host:</b>	Mouse
<b>Antigen:</b>	recombinant human TIF2
<b>Clone:</b>	3C11
<b>Isotype:</b>	IgG1, κ



<b>Application(s):</b>	ELISA
	Western Blot
	Immunofluorescence
	ChIP
	Recommended dilutions: 1/500-1/5000

**Specificity:** human TIF2. Epitope localized between aa 624 and 869 of human TIF2

### Scientific Literature

Duteil D, Chambon C, Ali F, Malivindi R, Zoll J, Kato S, Geny B, Chambon P, Metzger D. The transcriptional coregulators TIF2 and SRC-1 regulate energy homeostasis by modulating mitochondrial respiration in skeletal muscles. *Cell Metab.* 2010 Nov 3;12(5):49

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>