



## GCN5<sup>Δ</sup> Monoclonal Antibody

Cat No: G01045 - 100  $\mu$ L

### 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>	synthetic peptide coupled to ovalbumin
<b>Clone:</b>	2C11
<b>Isotype:</b>	IgG1, $\kappa$



NEW

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

**Specificity:** human and mouse GCN5-L

### Scientific Literature

Brand M, Moggs JG, Oulad-Abdelghani M, Lejeune F, Dilworth FJ, et al. (2001) UVdamaged DNA-binding protein in the TFIIIC complex links DNA damage recognition to nucleosome acetylation. *Embo J* 20: 3187-3196

Helmlinger D, Hardy S, Abou-Sleymane G, Eberlin A,

FP/25/20

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>