



## Gal-4 (DNA Binding domain) Monoclonal Antibody

Cat No: G01015 - 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>	Gal 4 DBD fusion protein
<b>Clone:</b>	GV10
<b>Isotype:</b>	IgG1, $\kappa$



<b>Application(s):</b>	ELISA
	Immunoblotting on recombinant proteins
	Immunocytochemistry on transfected cells
	Inhibits strongly DNA binding of Gal 4 DBD
	Transcription assays : see reference
	Immunoprecipitation : not tested
	recommended dilution: 1/1000-1/5000

**Specificity:** Gal 4 DBD containing proteins, epitope located in the first half of the DBD from aa 1 to 74

### Product Overview

This protein is a positive regulator for the gene expression of the galactose-induced genes such as GAL1, GAL2, GAL7, GAL10, and MEL1 which code for the enzymes used to convert galactose to glucose.

### Scientific Literature

White & al, EMBO J.(1992) vol.11, n°6 : 2229-2240

FP/19/20

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>