



Gal-4 (DNA Binding domain) Monoclonal Antibody

Cat No: G01016 - 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:	Gal 4 DBD fusion protein
Clone:	GV5
Isotype:	IgG1, κ



Application(s):	ELISA Immunoblotting on recombinant proteins Immunocytochemistry on transfected cells No supershift experiments Transcription assays : see reference Immunoprecipitation : not tested recommended dilution: 1/1000-1/5000
------------------------	---

Specificity: Gal 4 DBD fusion proteins, epitope located in the second half of the DBD from aa 74 to 147

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/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>