

Product Information

logo

ETS1 (DBD) Monoclonal Antibody

Cat No: G01014 - 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:	synthetic peptide coupled to ovalbumine
Clone:	7F2
Isotype:	IgG1, κ
Application(s):	ELISA Western blot Immunofluorescence recommended dilution: 1/500-1/5000
Specificity:	human and mouse ETS-1

Product Overview

ETS is one of the largest family of TF (with 28 members in Human). Protein C-ets-1 is a protein that in humans is encoded by the ETS1. ETS1 is playing an important role in immune cells, and tissues.

FP/18/20

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>