



MAT1-p32 / MNAT1 Monoclonal Antibody

Cat No: G01019 - 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:	Synthetic peptide
Clone:	1G6
Isotype:	IgG1, κ



Application(s):	ELISA
	Western Blot
	Immunocytochemistry
	recommended dilution: 1/500-1/5000

Specificity: human MAT1-p32.

Product Overview

MAT1 also known as MNAT1 belong to the complex of Cyclin dependent kinase-activating kinase, together with CDK-7 and Cyclin H where it plays the key role for regulating the activity of the complex. It is also playing a significant role in cell cycle, transcription and differentiation.

Scientific Literature

Busso D, Keriel A, Sandrock B, Poterszman A, Gileadi O, Egly JM. Distinct regions of MAT1 regulate cdk7 kinase and TFIIF transcription activities. J Biol Chem. 2000 Jul 28;275(30):22815-23.

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>