

# O(6)-Methylguanine-DNA Methyltransferase (MGMT) monoclonal antibody

This antibody stains all mantle zone lymphocytes and 50% of germinal center lymphocytes. Basaloid epithelial cells of tonsil squamous mucosa are also stained positive with this antibody.

MGMT, or O-6-methylguanine-DNA methyltransferase, is a crucial enzyme involved in DNA repair mechanisms within cells. Its primary function is to remove alkyl groups from the O-6 position of guanine, a common site of DNA damage induced by alkylating agents. By repairing this type of DNA damage, MGMT plays a vital role in maintaining genomic stability and protecting cells from the harmful effects of mutagenesis and carcinogenesis.

This antibody is covered by our [Worry-Free Guarantee](#).

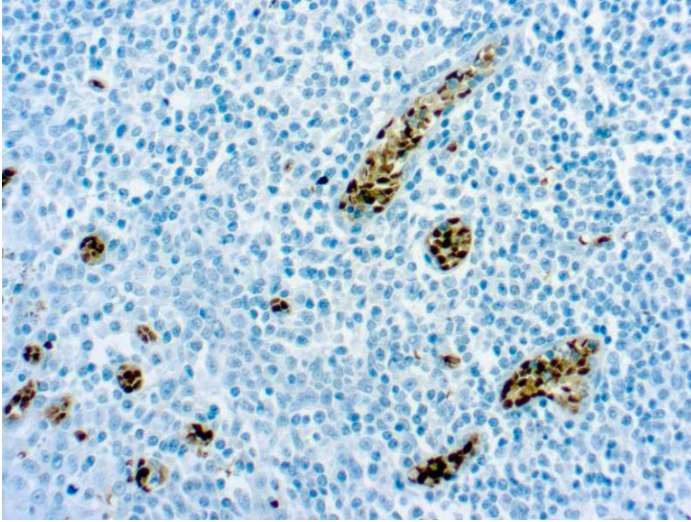
## Ordering Information

[Order Online »](#)

ENZ-ABS856-1000	1ml
ENZ-ABS856-0100	100µl

## Manuals, SDS & CofA

[View Online »](#)



Formalin fixed paraffin embedded human tonsil stained with MGMT antibody

## Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	MGMT
Application	IHC (PS)
Application Notes	Pretreatment: Citrate Buffer pH 6.0
Clone	MT3.1
Formulation	This product is supplied as an immunoglobulin and contains sodium azide as a preservative
Host	Mouse
Immunogen	Recombinant human MGMT protein
Isotype	IgG1
Positive Control	Tonsil
Recommendation Dilutions/Conditions	1:25-1:100Antibody dilutions should be appropriately adjusted and verified according to the detection system used.Optimal conditions must be determined individually for each application.
Species Reactivity	Human
Worry-free Guarantee	This antibody is covered by our <a href="#">Worry-Free Guarantee</a>