

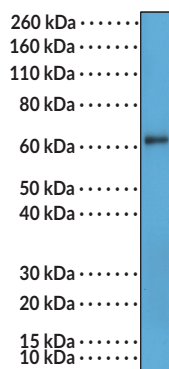
# PRODUCT INFORMATION

## Ecto-5'-nucleotidase/CD73 Rabbit Monoclonal Antibody (Clone RM431) Item No. 35613

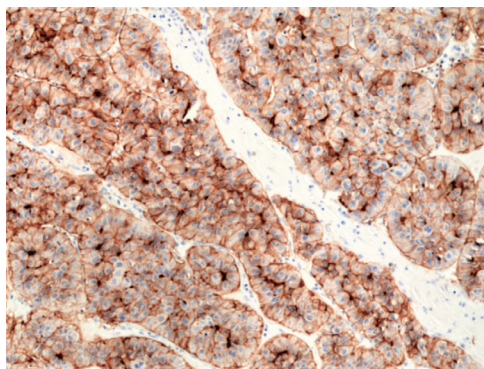
### Overview and Properties

<b>Contents:</b>	This vial contains 100 µl of protein A-affinity purified monoclonal antibody.
<b>Synonym:</b>	NT5E
<b>Immunogen:</b>	Recombinant human CD73
<b>Cross Reactivity:</b>	(+) CD73
<b>Species Reactivity:</b>	(+) Human
<b>Form:</b>	Liquid
<b>Storage:</b>	-20°C (as supplied)
<b>Stability:</b>	≥1 year
<b>Storage Buffer:</b>	PBS with 50% glycerol, 1% BSA, and 0.09% sodium azide
<b>Clone:</b>	RM431
<b>Host:</b>	Rabbit
<b>Isotype:</b>	IgG
<b>Applications:</b>	Immunohistochemistry (IHC) and Western blot (WB); the recommended starting dilution is 1:100-1:250 for IHC and 1:1,000-1:2,000 for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

### Images



WB of MDA-MB-231 cell lysate using CD73 Rabbit Monoclonal Antibody at a dilution of 1:1,000.



Immunohistochemical staining of formalin-fixed and paraffin-embedded human liver cancer tissue using CD73 Rabbit Monoclonal Antibody at a dilution of 1:200.

**WARNING**  
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

**SAFETY DATA**  
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

**WARRANTY AND LIMITATION OF REMEDY**  
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 01/17/2024

**CAYMAN CHEMICAL**  
1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA  
PHONE: [800] 364-9897  
[734] 971-3335  
FAX: [734] 971-3640  
CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM

# PRODUCT INFORMATION



## Description

---

CD73, also known as ecto-5'-nucleotidase, is a membrane-bound ectoenzyme that catalyzes the hydrolysis of AMP to adenosine, a mediator of purinergic signaling pathways.<sup>1,2</sup> It exists as a homodimer where each monomer is composed of a catalytic N-terminal domain, a C-terminal substrate-binding domain, and a glycosylphosphatidylinositol (GPI) anchor that is attached to the cell membrane.<sup>1</sup> CD73 is ubiquitously expressed, and a soluble form of CD73 (sCD73) can be produced by proteolytic cleavage of its GPI anchor.<sup>1</sup> It is upregulated by hypoxia and a variety of pro-inflammatory mediators, including TGF- $\beta$ , IFN, TNF- $\alpha$ , IL-1 $\beta$ , and prostaglandin E<sub>2</sub> (PGE<sub>2</sub>).<sup>3</sup> The enzymatic activity of CD73 is coupled to CD39, the enzyme that produces AMP from ATP, a danger signal produced from injured or dying cells.<sup>2</sup> CD73 converts AMP to adenosine, a ligand for adenosine receptors, which regulate various physiological responses, including inflammatory pathways.<sup>2,3</sup> It has other immunoregulatory functions, including acting as a T cell co-stimulatory molecule, inhibiting macrophage-mediated inflammation, and facilitating lymphocyte tethering to the endothelium.<sup>2</sup> CD73 has additional roles in the maintenance of tissue barrier function, cardioprotection during ischemia-reperfusion injury, and cancer, where it promotes metastasis, angiogenesis, and immune evasion.<sup>4</sup> Increased tumor CD73 expression is associated with poor overall survival in patients with breast or ovarian cancer.<sup>5</sup> Cayman's Ecto-5'-nucleotidase/CD73 Rabbit Monoclonal Antibody (Clone RM431) can be used for immunohistochemistry (IHC) and Western blot (WB) applications.

## References

---

1. Nocentini, A., Capasso, C., and Supuran, C.T. Small-molecule CD73 inhibitors for the immunotherapy of cancer: A patent and literature review (2017-present). *Expert Opin. Ther. Pat.* **31(10)**, 867-876 (2021).
2. Minor, M., Alcedo, K.P., Battaglia, R.A., et al. Cell type- and tissue-specific functions of ecto-5'-nucleotidase (CD73). *Am. J. Physiol. Cell Physiol.* **317(6)**, C1079-C1092 (2019).
3. Antonioli, L., Pacher, P., Vizi, E.S., et al. CD39 and CD73 in immunity and inflammation. *Trends Mol. Med.* **19(6)**, 355-367 (2013).
4. Gao, Z., Dong, K., and Zhang, H. The roles of CD73 in cancer. *Biomed. Res. Int.* 460654 (2014).
5. de Leve, S., Wirsdörfer, F., and Jendrossek, V. Targeting the immunomodulatory CD73/adenosine system to improve the therapeutic gain of radiotherapy. *Front. Immunol.* **10**, 698 (2019).

CAYMAN CHEMICAL  
1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA  
PHONE: [800] 364-9897  
[734] 971-3335  
FAX: [734] 971-3640  
CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM