

PRODUCT INFORMATION

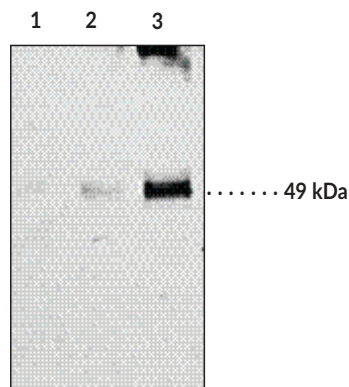


Adenosine Receptor A_{2A} Monoclonal Antibody (Clone 7F6-G₅-A2) Item No. 10011454

Overview and Properties

Contents:	This vial contains 500 µg of affinity-purified monoclonal antibody.
Synonyms:	A _{2A} R, ADORA2A
Immunogen:	Human full length A _{2A} R; the antibody recognizes amino acids 213-220 (SQPLPGER) as determined by epitope mapping
Cross Reactivity:	(+) Human A _{2A} R; other species not tested but the epitope is identical in rodents, primates, and canines and is expected to work for these samples. (-) A ₁ R, A _{2B} R
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥3 years
Storage Buffer:	PBS, pH 7.2, with 50% glycerol, 0.5 mg/ml BSA, and 0.02% sodium azide
Clone:	7F6-G ₅ -A2
Host:	Mouse
Isotype:	IgG2a
Applications:	Western blot (WB); the recommended starting concentration is 5 µg/ml. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Image



- Lane 1:** A_{2A}R transfected HEK cell membrane fraction (6.5 µg)
Lane 2: A_{2A}R transfected HEK cell membrane fraction (13 µg)
Lane 3: A_{2A}R transfected HEK cell membrane fraction (26 µg)

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, 11/01/2023

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

A_{2A}R is a multi-pass membrane protein that is normally localized to the plasma membrane.¹ This receptor is part of the G protein-coupled receptor family that binds adenosine and serves multiple functions. Antagonists of this receptor have been targeted for the treatment of Parkinson's disease.^{2,3} Early reports found this receptor is found primarily in the brain striatum, but is also found in immune cells and other tissues as well.² A_{2A}R is comprised of 412 amino acids with an expected molecular weight of 45 kDa.^{4,5} However multiple glycosylation sites exist that may explain the retarded migration observed by western blotting (45-50 kDa).¹ This antibody has been extensively characterized and the epitope has been mapped to the third intracellular loop of the receptor.⁵

References

1. Piersen, C.E., True, C.D., and Well, J.N. A carboxyl-terminally truncated mutant and nonglycosylated A_{2a} adenosine receptors retain ligand binding. *Mol. Pharmacol.* **45(5)**, 861-870 (1994).
2. Fredholm, B.B., Chern, Y., Franco, R., *et al.* Aspects of the general biology of adenosine A_{2A} signaling. *Prog. Neurobiol.* **83(5)**, 263-276 (2007).
3. Ledent, C., Vaugeois, J.-M., Schiffmann, S.N., *et al.* Aggressiveness, hypoalgesia and high blood pressure in mice lacking the adenosine A_{2a} receptor. *Nature* **388(6643)**, 674-678 (1997).
4. Robeva, A.S., Woodward, R.L., Jin, X., *et al.* Molecular characterization of recombinant human adenosine receptors. *Drug Development Research* **39**, 243-252 (1996).
5. Rosin, D.L., Robeva, A., Woodward, R.L., *et al.* Immunohistochemical localization of adenosine A_{2A} receptors in the rat central nervous system. *J. Comp. Neurol.* **401(2)**, 1631-186 (1998).

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