

# PRODUCT INFORMATION



## Adenosine A<sub>2A</sub> Receptor (human recombinant) Item No. 24730

### Overview and Properties

**Source:** N-terminal histidine- and strep-tagged human adenosine A<sub>2A</sub> receptor containing a TEV protease cleavage site purified from a baculovirus overexpression system in Sf9 cells

**Amino acids:** Full length, wild-type sequence

**Uniprot No.:** P29274

**Molecular Weight:** ~47.7 kDa

**Storage:** -80°C (as supplied)

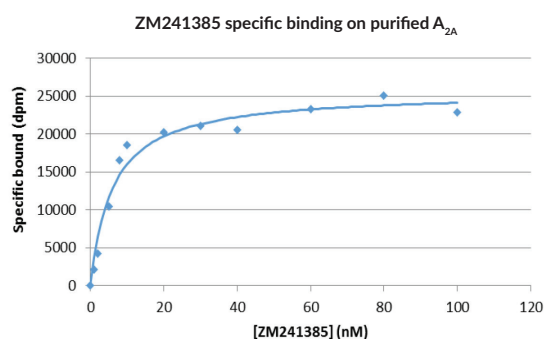
**Stability:** ≥2 years

**Purity:** *batch specific* (≥90% estimated by SDS-PAGE)

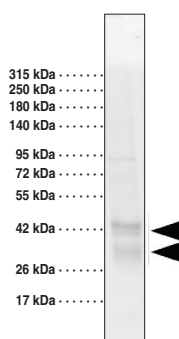
**Supplied in:** 50 mM HEPES, pH 7.4, 200 mM sodium chloride, and 0.05%/0.006% DDM/CHS

*Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.*

### Images



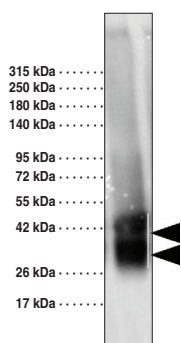
QC: Activity measured by radiobinding assay. Binding of [<sup>3</sup>H]ZM241385 was measured on purified A<sub>2A</sub> receptor. A K<sub>d</sub> of 6 nM was determined for ZM241385.



SDS-PAGE, 4-15% acrylamide gel Bio-rad Stain-Free™ detection

SDS-PAGE, Stain-Free™ detection

Purified A<sub>2A</sub> receptor was migrated on a 4-15% Tris-glycine SDS-PAGE and the total proteins were Stain-Free™ detected. Black arrows indicate the target. Upper arrow indicates full length A<sub>2A</sub> receptor. Lower arrow indicates shorter A<sub>2A</sub> receptor resulting from partial cleavage of C-term end.



SDS-PAGE, 4-15% acrylamide gel WB Anti-ICL3 A<sub>2A</sub> antibody (7F6-G5-A2)

SDS-PAGE, Western blotting

Purified A<sub>2A</sub> receptor was migrated on a 4-15% Tris-glycine SDS-PAGE, transferred to PVDF membrane and immunodetected with a monoclonal anti-ICL3-A<sub>2A</sub> (7F6-G5-A2, SCB7). Black arrows indicate the target.

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, 05/25/2018

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

---

Adenosine A<sub>2A</sub> receptor (human recombinant) contains amino acids corresponding to the full length, wild-type, human A<sub>2A</sub> receptor, as well as N-terminal histidine and streptavidin tags with a TEV protease cleavage site. The A<sub>2A</sub> receptor is a G protein-coupled receptor (GPCR) that is activated by adenosine on presynaptic terminals to induce neurotransmitter release and on postsynaptic terminals to increase excitability.<sup>1</sup> Signaling through the A<sub>2A</sub> receptor occurs through G<sub>s</sub>, G<sub>olf</sub>, and G<sub>15/16</sub> proteins to increase cyclic AMP or inositol-(1,4,5)-trisphosphate (IP<sub>3</sub>) activity.<sup>2</sup> It is expressed in the brain on striatal-pallidal GABAergic neurons and in the olfactory bulb as well as other brain regions. It is also expressed in the periphery in organs and cells involved with immune function, including the spleen, thymus, leukocytes, and blood platelets as well as in the heart, lung, and blood vessels. A<sub>2A</sub> receptor activity induced by agonists has neurodegenerative effects, which has led to the development of A<sub>2A</sub> receptor antagonists, but has shown mixed effects on seizure activity in rodent models.<sup>1</sup> The A<sub>2A</sub> receptor also plays a role in the regulation of enteric nervous system function, vasodilation of coronary arteries, and the immune response.<sup>3</sup> In addition to signaling through the A<sub>2A</sub> receptor alone, it can form heteromers with other GPCRs, including adenosine A<sub>1</sub>, dopamine D<sub>2</sub> and D<sub>3</sub>, cannabinoid CB<sub>1</sub>, and metabotropic glutamate receptor 5 (mGluR5).<sup>4,5</sup> The function of these receptors is influenced by the interaction, with the A<sub>2A</sub> receptor decreasing dopamine binding to the D<sub>2</sub> receptor and mGluR5 potentiating the effects of the A<sub>2A</sub> receptor on adenylyl cyclase and MAPK signaling.<sup>4</sup>

## References

---

1. Stockwell, J., Jakova, E., and Cayabyab, F.S. Adenosine A1 and A2A receptors in the brain: Current research and their role in neurodegeneration. *Molecules* **22**(4), pii: E676 (2017).
2. Fredholm, B.B., Arslan, G., Halldner, L., et al. Structure and function of adenosine receptors and their genes. *Naunyn Schmiedebergs Arch. Pharmacol.* **362**(4-5), 364-374 (2000).
3. Sousa, J.B. and Diniz, C. The adenosinergic system as a therapeutic target in the vasculature: New ligands and challenges. *Molecules* **22**(5), pii: E752 (2017).
4. Ferré, S., Ciruela, F., Quiroz, C., et al. Adenosine receptor heteromers and their integrative role in striatal function. *ScientificWorldJournal* **7**, 74-85 (2007).
5. Torvinen, M., Marcellino, D., Canals, M., et al. Adenosine A<sub>2A</sub> receptor and dopamine D<sub>3</sub> receptor interactions: Evidence of functional A<sub>2A</sub>/D<sub>3</sub> heteromeric complexes. *Mol. Pharmacol.* **67**(2), 400-407 (2005).

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