

# Nucleoside and Nucleotide Ligands for Nucleoside Receptors (A<sub>1</sub>, A<sub>2A</sub> and A<sub>3</sub>)

Nucleosides and Nucleotides	Cat. No.	Subreceptor Type
<b>Recommended Ligands for Subreceptor Type A<sub>1</sub></b>		
<b>N<sup>6</sup>-Cyclopentyl-adenosine (CPA)</b>	N-1078	A <sub>1</sub> , A <sub>2A</sub> , A <sub>3</sub>
<b>5'-Ethylcarboxamido-adenosine (NECA)</b>	N-1076	A <sub>1</sub> , A <sub>2A</sub> , A <sub>3</sub>
<b>N<sup>6</sup>-Methyl-adenosine</b>	N-1080	A <sub>1</sub> , A <sub>2A</sub> , A <sub>3</sub>
<b>2-Chloro-N<sup>6</sup>-cyclopentyl-adenosine (CCPA)</b>	N-1079	A <sub>1</sub>
<b>8-Cyclopentyl-1,3-dipropyl-xanthine (DPCPX)</b>	N-1077	A <sub>1</sub>
<b>Recommended Ligands for Subreceptor Type A<sub>2A</sub></b>		
<b>N<sup>6</sup>-Cyclopentyl-adenosine (CPA)</b>	N-1078	A <sub>1</sub> , A <sub>2A</sub> , A <sub>3</sub>
<b>5'-Ethylcarboxamido-adenosine (NECA)</b>	N-1076	A <sub>1</sub> , A <sub>2A</sub> , A <sub>3</sub>
<b>N<sup>6</sup>-Methyl-adenosine</b>	N-1080	A <sub>1</sub> , A <sub>2A</sub> , A <sub>3</sub>
<b>Recommended Ligands for Subreceptor Type A<sub>3</sub></b>		
<b>N<sup>6</sup>-Cyclopentyl-adenosine (CPA)</b>	N-1078	A <sub>1</sub> , A <sub>2A</sub> , A <sub>3</sub>
<b>5'-Ethylcarboxamido-adenosine (NECA)</b>	N-1076	A <sub>1</sub> , A <sub>2A</sub> , A <sub>3</sub>
<b>N<sup>6</sup>-Methyl-adenosine</b>	N-1080	A <sub>1</sub> , A <sub>2A</sub> , A <sub>3</sub>
<b>N<sup>6</sup>-(2-Phenylethyl)-adenosine</b>	N-1080	A <sub>3</sub>
<b>Chloro-IB-MECA</b>	N-1075	A <sub>3</sub>
<b>Analogues of Adenosines</b>		
<b>β-L-Adenosine</b>	N-RP-9031	A <sub>1</sub> , A <sub>3</sub>
<b>N<sup>6</sup>-Benzylaminopurine riboside</b>	N-1022	A <sub>1</sub>
<b>8-Bromo-adenosine</b>	N-RP-1214	A <sub>1</sub>
<b>Arabinofuranosyl-adenine (ara-A)</b>	N-20305	A <sub>1</sub> , A <sub>2A</sub> , A <sub>3</sub>
<b>2'-Deoxyadenosine monohydrate</b>	N-DN-1001	A <sub>3</sub>
<b>Analogues of ATP</b>		
<b>ATP</b>	NU-1010	A <sub>1</sub>
<b>NPE-caged-ATP</b>	NU-301	A <sub>1</sub> release technology
<b>DMB-caged-ATP</b>	NU-309	A <sub>1</sub> release technology
<b>ATPyS</b>	NU-406	A <sub>1</sub>
<b>AppCp</b>	NU-422	A <sub>1</sub>