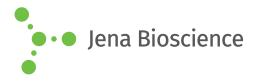
DATA SHEET

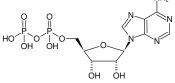




ADP - Solid

Adenosine-5'-diphosphate, Sodium salt

Cat. No.	Amount
NU-1198-1G	1 g
NU-1198-5G	5 g



For general laboratory use.

Shipping: shipped on gel packs

Storage Conditions: store at -20 °C

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

Shelf Life: 12 months after date of delivery

Molecular Formula: $C_{10}H_{15}N_5O_{10}P_2$ (free acid)

Molecular Weight: 427.20 g/mol (free acid)

Exact Mass: 427.03 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: solid

Color: white to off-white

Spectroscopic Properties: λ_{max} 259 nm, ϵ 15.4 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

Applications:

Agonistic ligand, mainly for nucleoside receptor A1

Nucleoside-triphosphates can be converted by different membranebound phosphatases into nucleosides acting as nucleoside receptor ligands. In some cases nucleoside phosphates act also directly on nucleoside receptors.

Selected References:

Volonte *et al.* (2009) Membrane components and purinergic signalling: the purinome, a complex interplay among ligands, degrading enzymes, receptors and transporters. *FEBS J.* **276**:318.

Yegutkin (2008) Nucleotide and nucleoside converting enzymes: Important modulators of purinergic signalling cascade. *Biochim. Biophys. Acta* **1783**:673.

Williams *et al.* (1986) Effects of purine nucleotides on the binding of [3H]cyclopentyladenosine to adenosine A1-receptors in rat brain membranes. *J. Neurochem.* **47 (1)**:88.

