

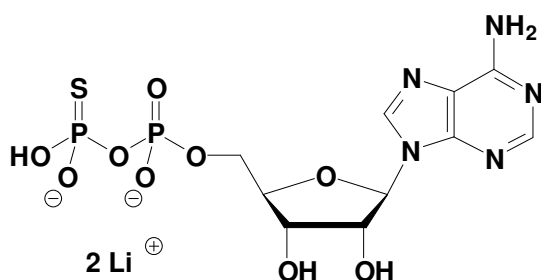
## ADP $\beta$ S

Adenosine-5'-( $\beta$ -thio)-diphosphate, Lithium salt

Cat. No.	Amount
NU-433S	10 mg
NU-433L	25 mg

### Selected References:

Goody *et al.* (1971) Thiophosphate Analogs of Nucleoside Di- and Triphosphates. *J. Amer. Chem. Soc.* **93**:6252.



**Cat. No.:** NU-433

**Molecular Formula:** C<sub>10</sub>H<sub>13</sub>N<sub>5</sub>O<sub>9</sub>P<sub>2</sub>S (Anion)

**Molecular Weight:** 441.24 (Anion)

**Purity:** > 85%, lyophilized powder

### Storage conditions:

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at < -20°C. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 6 months after date of delivery.

**For research use only!**