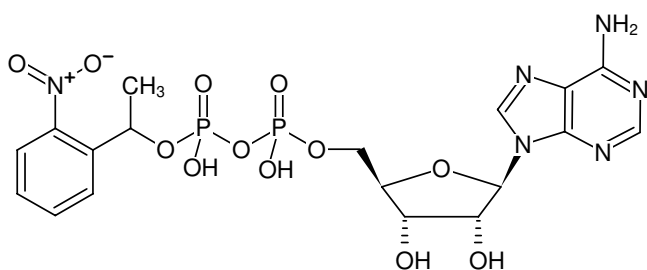


**NPE-caged-ADP**Adenosine-5'-diphosphate, P<sup>2</sup>-(1-(2-nitrophenyl)-ethyl)-ester, Potassium salt

Cat. No.	Amount
NU-952S	150 µl (10 mM)
NU-952L	5 x 150 µl (10 mM)



Structural formula of NPE-caged-ADP

**For general laboratory use.****Shipping:** shipped on gel packs**Storage Conditions:** store at -20 °C**Additional Storage Conditions:** store dark

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

**Shelf Life:** 12 months after date of delivery**Molecular Formula:** C<sub>18</sub>H<sub>22</sub>N<sub>6</sub>O<sub>12</sub>P<sub>2</sub> (free acid)**Molecular Weight:** 576.35 g/mol (free acid)**Exact Mass:** 576.077 g/mol (free acid)**CAS#:** 340830-06-0 (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** solution in water**Color:** colorless to slightly yellow**Concentration:** 10 mM - 11 mM**pH:** 7.5 ±0.5**Spectroscopic Properties:** λ<sub>max</sub> 260 nm, ε 18.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)