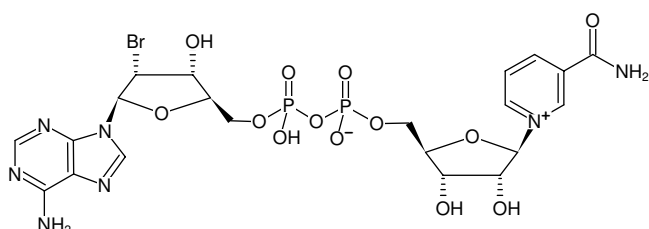


**2'-Bromo-dAP<sub>2</sub>(Nic)**(2'Br-dAP<sub>2</sub>(Nic))P<sup>1</sup>-[5'-(2'-Bromo-2'-deoxy-adenosyl)] P<sup>2</sup>-[5'-(1-β D-Ribofuranosyl-nicotinamide)] diphosphate, Sodium salt

Cat. No.	Amount
NU-524	300 µl (10 mM)

Structural formula of 2'-Bromo-dAP<sub>2</sub>(Nic)**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<sub>21</sub>H<sub>26</sub>BrN<sub>7</sub>O<sub>13</sub>P<sub>2</sub> (free acid)**Molecular Weight:** 726.32 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** solution in water**Color:** colorless to slightly yellow**Concentration:** 10 mM - 11 mM**pH:** 7.5 ±0.5