

V. 1.0

Adenosine 5'-triphosphate disodium salt (ATP)

Cat. # BSA04

Upon arrival store at 4°C (desiccated) See datasheet for storage after reconstitution

Material

Adenosine 5'-triphosphate (ATP) is supplied as a white powder. The compound is derived from a microbial source. The molecular weight of the compound is 551.14.

Chemical formula $C_{10}H_{14}N_5Na_2O_{13}P_3$. xH_2O . CAS # 42373-41-1.

Storage and Reconstitution

Briefly centrifuge to collect the product at the bottom of the tube. The lyophilized compound, when stored desiccated to <10% humidity at 4°C is stable for 6 months. The ATP should be reconstituted to 100 mM with 1 ml of cold 100mM Tris pH 7.5. The ATP should then be aliquoted into experiment sized amounts, snap frozen in liquid nitrogen and stored at or below -20°C. The ATP stock is stable for 6 months if stored at or below -20°C.

Purity

Compound is greater than 99% pure as determined by HPLC analysis.

Product Citations / Related Products

For the latest citations and related products please visit www.cytoskeleton.com.