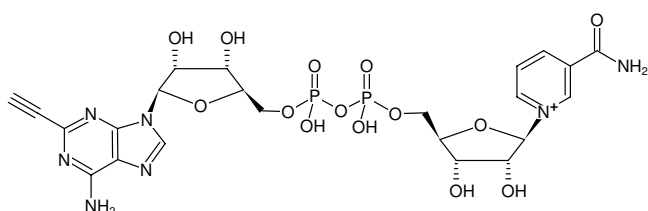




## 2-Ethynyl-Adenosine-NAD<sup>+</sup>

(2-EA-NAD<sup>+</sup>)

Cat. No.	Amount
CLK-043	20 µl (10 mM)



Structural formula of 2-Ethynyl-Adenosine-NAD<sup>+</sup>

### For research use only!

**Shipping:** shipped on dry ice

**Storage Conditions:** store at -20 °C

**Shelf Life:** 12 months after date of delivery

**Molecular Formula:** C<sub>23</sub>H<sub>28</sub>N<sub>7</sub>O<sub>14</sub>P<sub>2</sub> (cation)

**Molecular Weight:** 688.46 g/mol (cation)

**Exact Mass:** 688.12 g/mol (cation)

**Form:** solution in water

**Color:** colorless to slightly yellow

**Concentration:** 10 mM - 11 mM

### Applications:

*in vitro* ADP-Ribosylation of proteins<sup>[1](compound 1)</sup>

The resulting alkyne-functionalized protein<sup>[1,2]</sup> can subsequently be processed via Cu(I)-catalyzed (azide-alkyne) click chemistry that offers the choice

- to introduce a Biotin group for subsequent purification tasks (via Azides of Biotin)
- to introduce fluorescent group for subsequent microscopic imaging (via Azides of fluorescent dyes)

Presolski *et al.*<sup>[3]</sup> and Hong *et al.*<sup>[4]</sup> provide a general protocol for Cu(I)-catalyzed click chemistry reactions and Lin *et al.*<sup>[2]</sup> for ADP-Ribosylation assays that may be used as a starting point for the set up and optimization of individual assays.

### Related Products:

Copper (II)-Sulphate (CuSO<sub>4</sub>), #CLK-MI004

Tris(3-hydroxypropyltriazolylmethyl)amine (THPTA), #CLK-1010

Sodium Ascorbate (Na-Ascorbate), #CLK-MI005

### Selected References:

[1] Wang *et al.* (2014) Chain-Terminating and Clickable NAD<sup>+</sup> analogues for labeling the target proteins of ADP-Ribosyltransferases. *Angew. Chem. Int. Ed.* **53**:8159.

[2] Lin *et al.* (2012) Labeling substrate proteins of Poly (ADP-ribose) polymerase with Clickable NAD analog. *Curr. Prot. in Chem. Biol.* **4**:19.

[3] Presolski *et al.* (2011) Copper-Catalyzed Azide-Alkyne Click Chemistry for Bioconjugation. *Current Protocols in Chemical Biology* **3**:153.

[4] Hong *et al.* (2011) Analysis and Optimization of Copper-Catalyzed Azide-Alkyne Cycloaddition for Bioconjugation. *Angew. Chem. Int. Ed.* **48**:9879.