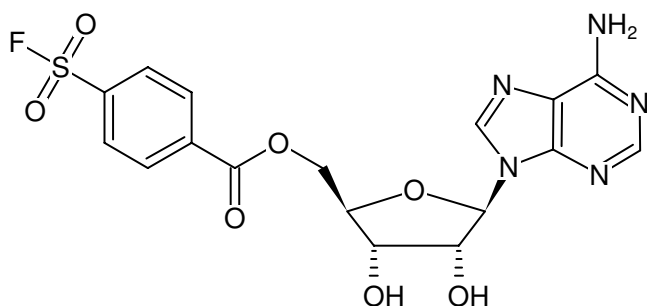


**5'-O-(4-fluorosulfonylbenzoyl)-adenosine**

(FSBA)

Cat. No.	Amount
N-1086-25	25 mg



Structural formula of 5'-O-(4-fluorosulfonylbenzoyl)-adenosine

**For general laboratory use.****Shipping:** shipped at ambient temperature**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>17</sub>H<sub>16</sub>FN<sub>5</sub>O<sub>7</sub>S**Molecular Weight:** 453.40 g/mol**Exact Mass:** 453.08 g/mol**Purity:** ≥ 95 % (HPLC)**Form:** solid**Color:** white to off-white**Selected References:**Embogama *et al.* (2017) K-BILDS: A Kinase Substrate Discovery Tool. *Chembiochem.* **18** (1):136.Knight *et al.* (2013) Profiling the kinome: current capabilities and future challenges. *J. Proteomics.* **81**:43.Knight *et al.* (2012) A novel whole-cell lysate kinase assay identifies substrates of the p38 MAPK in differentiating myoblasts *Skelet Muscle.*

doi:10.1186/2044-5040-2-5.