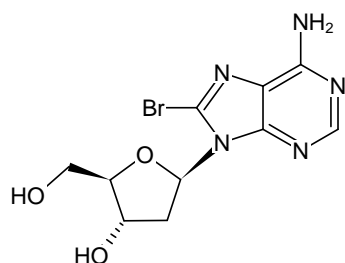


**8-Bromo-2'-deoxyadenosine**

Cat. No.	Amount
N-DN-1182-1G	1 g



Structural formula of 8-Bromo-2'-deoxyadenosine

***For research use only!*****Shipping:** shipped at ambient temperature**Storage Conditions:** store at -20 °C**Shelf Life:** 30 months after date of delivery**Molecular Formula:** C<sub>10</sub>H<sub>12</sub>BrN<sub>5</sub>O<sub>3</sub>**Molecular Weight:** 330.14 g/mol**Purity:** ≥ 98 % (HPLC)**Form:** white to off-white powder**Solubility:** 0.1 N HCl