

8-Br-Ade

8-Br-Ade can be used as a reactive precursor for the synthesis of 8-modified adenine analogues.

BLG-B052-50 (5 x 10 μ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

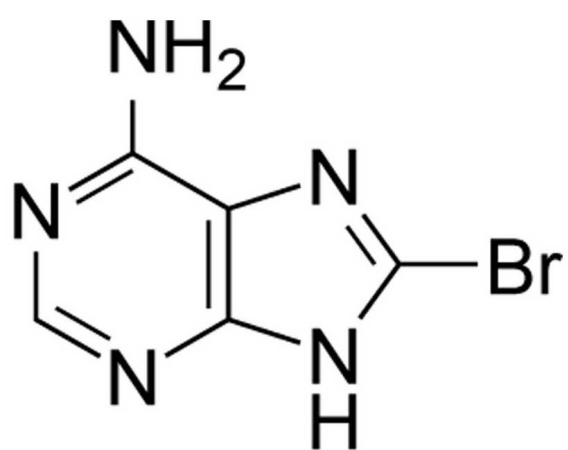
Ordering Information

[Order Online »](#)

| | |
|-------------|----------------|
| BLG-B052-50 | 5x10 μ mol |
|-------------|----------------|

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 8-Bromoadenine

CAS 6974-78-3

Couple Type Modified nucleotides

Formula $C_5H_4BrN_5$

MW 214

Purity > 95% HPLC

Quantity 10µmol≈2.1mg

Solubility Soluble in DMSO or DMF.

Technical Info / Product Notes For the Original Manufacturer's data sheet please [click here](#).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com