

5-Hydroxymethyl-dC (100 mg)

Ordering Information

[Order Online »](#)

JBS-N-1070-100

100mg

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 hmdC, 5-Hydroxymethyl-2'-deoxycytidine

Appearance white to off-white solid

Application Notes Epigenetic therapy
Metabolism of dC substituted at position 5
Influence on DNA promotor

CAS 7226-77-9

Formula $C_{10}H_{15}N_3O_5$

MW 257.24

Purity ≥ 97 % (HPLC)

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