
5-Azacytidine

(Catalog No. M22001)

Alternative Names

Azacytidine Vidaza

Description

DNA methyltransferase inhibitor which results in DNA demethylation (hypomethylation) and gene activation by remodeling "opening" chromatin. Genes are synergistically reactivated when demethylation is combined with histone hyperacetylation.

CAS Number

320-67-2

Chemical Formula

$C_8H_{12}N_4O_5$

Molecular Weight

244.2

Size

1 g

Purity

min 98%

Solubility

50% Acetic acid or DMSO

Storage Temp

-20°C

Ordering Information:

Products	Size	Cat. No.
5-Azacytidine	1 g	M22001-1

This product is for research purposes only. Not intended for use in diagnostic procedures.