

5-Methyl-CTP NU-1138-THR

Product Name

Long Name
Catalog Number
LOT Number
Formula
Molecular Weight
Exact Mass
Spectroscopic Properties
Storage
Shelf Life

5-Methyl-CTP

5-Methylcytidine-5'-triphosphate, Sodium Salt
NU-1138-THR
XXX
 $C_{10}H_{18}N_3O_{14}P_3$ (free acid)
 $497.18 \text{ g x mol}^{-1}$ (free acid)
 $497.00 \text{ g x mol}^{-1}$ (free acid)
 $\lambda_{\text{max}} = 277 \text{ nm}$; $\epsilon = 9.0 \text{ L x mmol}^{-1} \text{ x cm}^{-1}$ (Tris-HCl, pH 7.5)
store at $-20 \text{ }^\circ\text{C}$
12 months after date of delivery

Test

Form
Color
Concentration
pH
Purity
Identification
Functionality (T7 IVT, 1400 nt transcript, fluorescence microplate assay)
DNAses (gel electrophoresis)
RNAses (gel electrophoresis)
Protease (UV-Vis)
Endotoxin (LAL test)

Specification

solution in water
colorless to slightly yellow
100 mM – 105 mM
 7.5 ± 0.25
 $\geq 99 \%$ (RP-HPLC)
MS negative and/or positive mode
 $\geq 100 \mu\text{g RNA}$ (100% substitution)
not detectable (LOD: 0.002 U DNase I equiv.)
not detectable (LOD: 1 pg RNase A equiv.)
not detectable (LOD: 5 ng Proteinase K equiv.)
 $\leq 1.0 \text{ EU/ml}$

Result

conforms
conforms
conforms
conforms
XY %
conforms
conforms
conforms
conforms
conforms