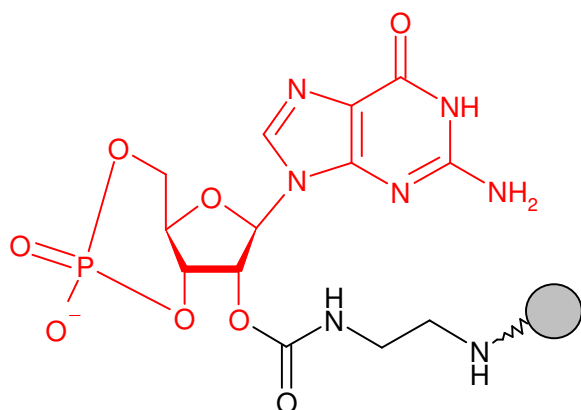


2'-EDA-cGMP - Sepharose®

Cat. No.	Amount
AC-148S	1 ml (bulk material)
AC-148L	5 ml (bulk material)
AC-148C	1 ml (syringe column)
AC-148SC	0.2 ml (scr. column)
AC-148MP	1 ml (MPLC column)



2'-O-(2-Aminoethyl-carbamoyl)-guanosine
-3',5'-cyclic-monophosphate, EDA-cGMP immobilized on
Agarose,

Degree of substitution: 5 μ mol EDA-cGMP/ml Gel 5
5 μ mol 2'/3'-EDA-m7-GDP/ml Gel

Linker: none

Storage buffer: 20% Ethanol

Storage: Short term exposure (up to 1 week cumulative) to ambient temperature possible.

Long term storage at 4°C.

Expiry: 12 month

® Sepharose is a trademark of Amersham-Biosciences.