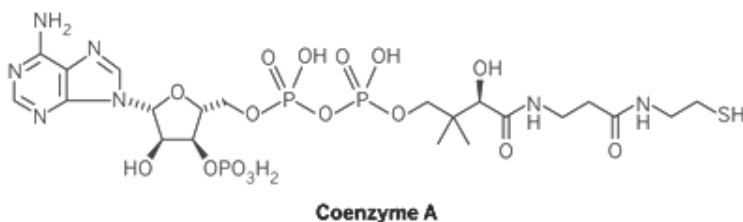


## Certificate of Analysis

### CAT# 309 Coenzyme A

Date: Mar 30, 2019  
 Compound ID: Coenzyme A (SH-CoA or CoA-SH)  
 CAS#: 85-61-0  
 Compound Name: Adenosyl-3-phospho-5-diphospho-pantothenate (reduced)  
 Batch Number: 5587  
 Molecular Formula: C<sub>21</sub>H<sub>36</sub>N<sub>7</sub>O<sub>16</sub>P<sub>3</sub>S  
 MW: 767.5 g/mol

Structure:



Test Items	Specification	Results	Test method
<b>Appearance</b>	Yellowish Solid	Conforms	Visual
<b>Loss on Drying</b>	<5%	1.3%	Weight
<b>Identification</b>	Activity consistent with structure	278 U/mg	> 250 U/mg
<b>Storage</b>	Store dry and dark at -20°C or for long term storage at -80°C		