

Certificate of Analysis

CAT#306 D-Luciferin

Date of Test: January, 17 2019

Compound ID: D-Luciferin (Firefly Luciferin)

CAS#: 2591-17-5

IUPAC Name: (4S)-2-(6-hydroxy-1,3-benzothiazol-2-yl)-4,5-dihydrothiazole-4-carboxylic acid

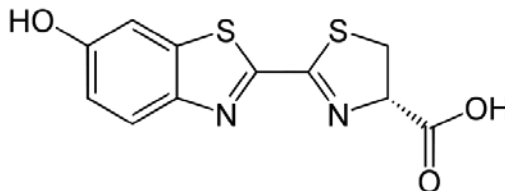
Batch Number: 5425

Manufacturer: Prolume Ltd.

Molecular Formula: C₁₁H₈N₂O₃S₂

MW: 280.32 g/mol

Structure:



Test Items	Specification	Results	Test method
Appearance	Light Yellow Solid	Conforms	Visual
Purity	>99%	99.091%	LCMS
Identification	Activity consistent with structure	Conforms	Light emission
	Spectrum is consistent with structure	Conforms	By HNMR
Storage	Store dry and dark at -20°C or for long term storage at -80°C.		