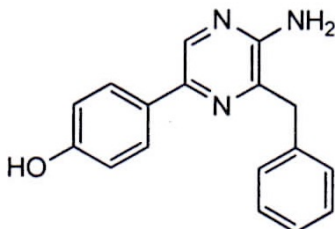


## Certificate of Analysis

### CAT #300 Coelenteramine

Date of Test: 22 May 2019  
 Compound ID: Coelenteramine (CLN)  
 CAS#: 37156-84-6  
 IUPAC Name: 4-(5-amino-6-benzylpyrazin-2-yl)phenol  
 Batch Number: 11711-012  
 Manufacturer: Prolume Ltd  
 Molecular Formula: C<sub>17</sub>H<sub>15</sub>N<sub>3</sub>O  
 MW: 277.32 g/mol

Structure:



Test Items	Specification	Results	Test method
Appearance	Yellow Solid	Conforms	Visual
Purity	>95%	96.98%	HPLC
Identification	Spectrum is consistent with structure	Conforms	By NMR
	Spectrum is consistent with structure	Conforms	By HPLC
Storage	Store at -20°C or for long term storage at -80°C.		