



3050 Spruce Street  
Saint Louis, Missouri 63103 USA  
Telephone 800-325-5832 • (314) 771-5765  
Fax (314) 286-7828  
email: techserv@sia.com  
sigma-aldrich.com

## Product Information

### COCAINE HYDROCHLORIDE

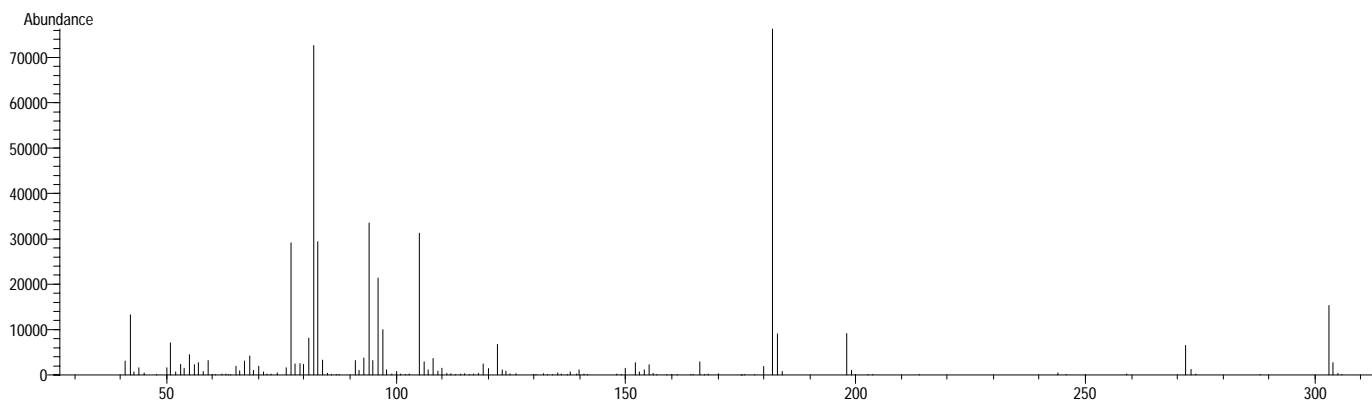
Methanol Solution

Product Number **C1528**

### TECHNICAL BULLETIN

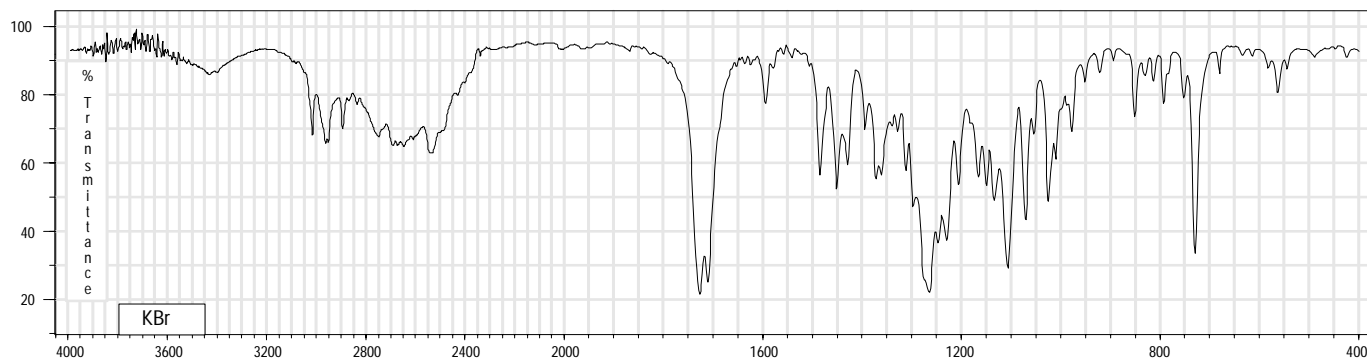
<b>Lot Number</b>	93K8800	<b>Synonyms</b>	3-(Benzoyloxy)-8-azabicyclo-[3.2.1]octane-2-carboxylic acid methyl ester hydrochloride;
<b>CAS Number</b>	[53-21-4]		l-Cocaine hydrochloride; Ecgonine
<b>Concentration</b>	1.12 mg Cocaine Hydrochloride (equivalent to 1.00 mg free base) /mL GC grade methanol (conc. verified by UV)		methyl ester bezoate hydrochloride; Benzoylmethylecgonine hydrochloride
<b>Molecular Formula</b>	C <sub>17</sub> H <sub>21</sub> NO <sub>4</sub> HCl	<b>Stability</b>	If stored at 2-8 °C in an air tight container, protected from light, decomposition of the product is less than 1% in 24 months. After opening, concentration may change due to loss of solvent.
<b>Molecular Weight</b>	339.8		
<b>DEA Class</b>	Exempt preparation of Class II		
<b>Expiration Date</b>	June 2006		
<b>Storage</b>	2-8 °C; After opening,		

70 eV Electron Ionization Mass Spectrum



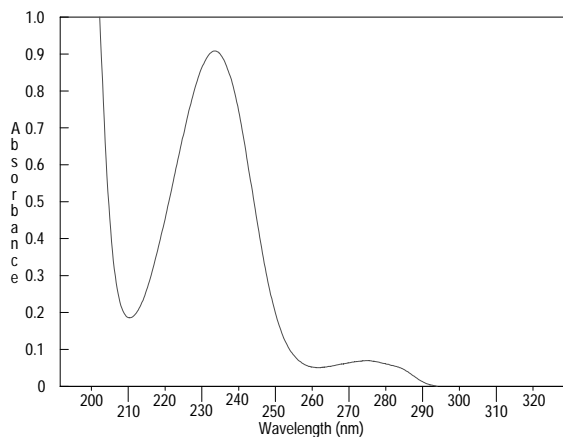
Hewlett-Packard 5890 Series II Gas Chromatograph  
Hewlett-Packard 5971 Mass Selective Detector  
Hewlett-Packard G1034C MS ChemStation (HP-DOS series)

## FT-IR



## Ultraviolet Spectrum

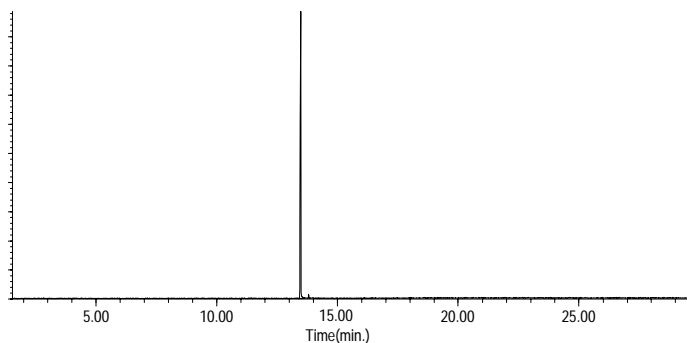
<b>Peak</b>	233.3 nm
<b>Absorbance</b>	0.908
<b>Solvent</b>	0.2 N H <sub>2</sub> SO <sub>4</sub>
<b>Dilution</b>	1 : 49 (v/v)



Bio-Tek Uvikon XL

## Capillary Gas Chromatography

**Column** Supelco SPB-1 (Cat. No. 2-4162)  
**I.D.** 0.20 mm **Length** 15 mm **Film Thickness** 0.20 µm  
**Oven Temperature** 50 °C (2 min.), then 15°/min. to 300 °C, hold 11 min.  
**Injector Temperature** 250 °C, Split  
**Detector** Hewlett-Packard 5971  
Mass Selective Detector (70 eV, 40-500 m/z)



Hewlett-Packard 5890

This product is not for *in vitro* diagnostic procedures. For R&D use only. Not for drug, household or other uses.  
Sigma brand products are sold through Sigma-Aldrich, Inc.  
Sigma-Aldrich, Inc. warrants that its products conform to the information contained in this and other Sigma-Aldrich publications.  
Purchaser must determine the suitability of the product(s) for their particular use. Additional terms and conditions may apply.  
Please see reverse side of the invoice or packing slip.