



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

## Product Information

### $\Delta^9$ -Tetrahydrocannabinol

Methanol Solution

**Product Number** T 4764  
**Lot Number** 93K8811  
**CAS Number** [1972-08-3]  
**Concentration** 1.03 mg  $\Delta^9$ -  
Tetrahydrocannabinol/mL GC  
grade methanol (conc. verified  
by HPLC)  
**Molecular Formula**  $C_{21}H_{30}O_2$   
**Molecular Weight** 314.5  
**DEA Class** Exempt  
preparation of

**Synonyms** Tetrahydro-669-trimethyl-3-pentyl-  
6H-dibenzo[bd]pyran-1-ol;  
 $\Delta^1$ -Tetrahydrocannabinol;  $\Delta^9$ -THC;  
 $\Delta^1$ -THC

**Stability** 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.

**Expiration Date** May 2006

#### High Pressure Liquid Chromatography

**Storage** 2-8 °C; After opening,  
transfer contents to small vial provided

**Column** C-18: 5 micron, 4.6 x 25 cm (Supelco 58298-C46)

**Mobile Phase**

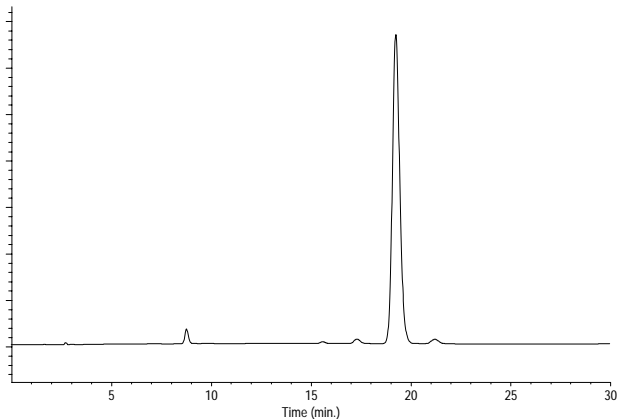
80 % Methanol / 20 % 18 M $\Omega$  water

**Gradient** isocratic

**Flow Rate** 1.0 mL/min.

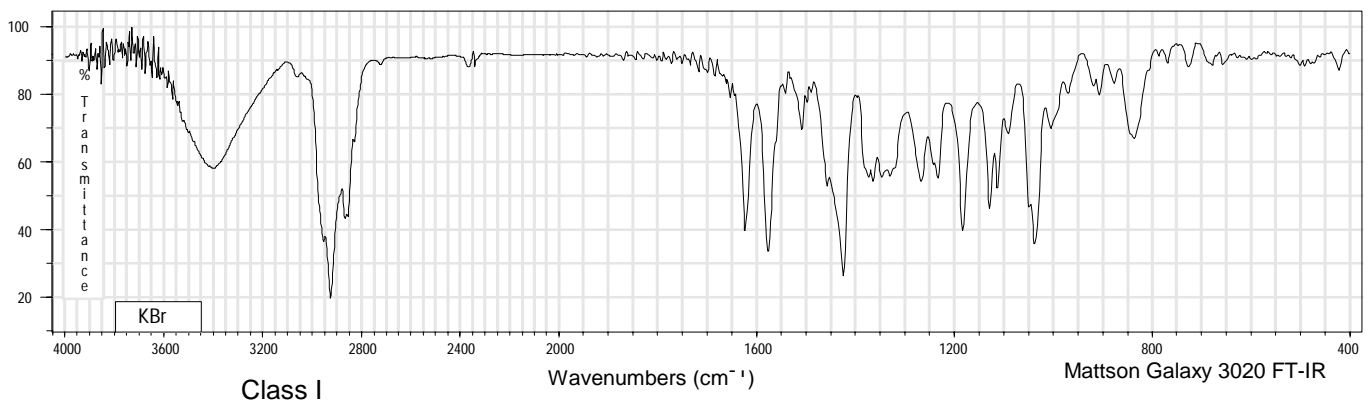
**Temp** ambient

**Detector** UV @ 210 nm



Perkin-Elmer LC-4 with Hewlett-Packard 3365 Series II Chemstation  
for your convenience.

#### FT-IR

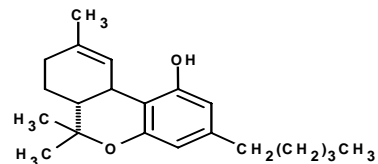


# $\Delta^9$ -TETRAHYDROCANNABINOL

Methanol Solution

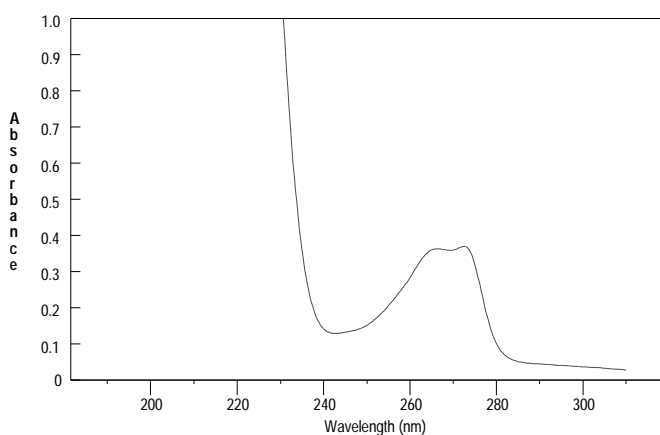
Product Number T 4764

Lot Number 93K8811



## Ultraviolet Spectrum

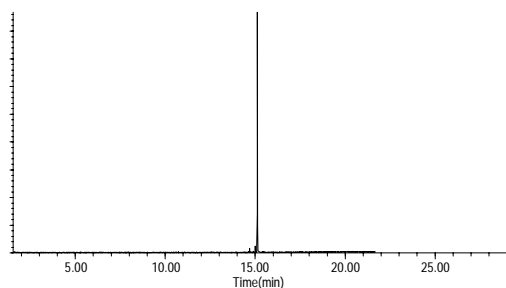
**Peak** 276.3nm  
**Absorbance** 0.363  
**Solvent** Methanol  
**Dilution** 1 : 9 (v/v)



Bio-Tek Uvikon XL

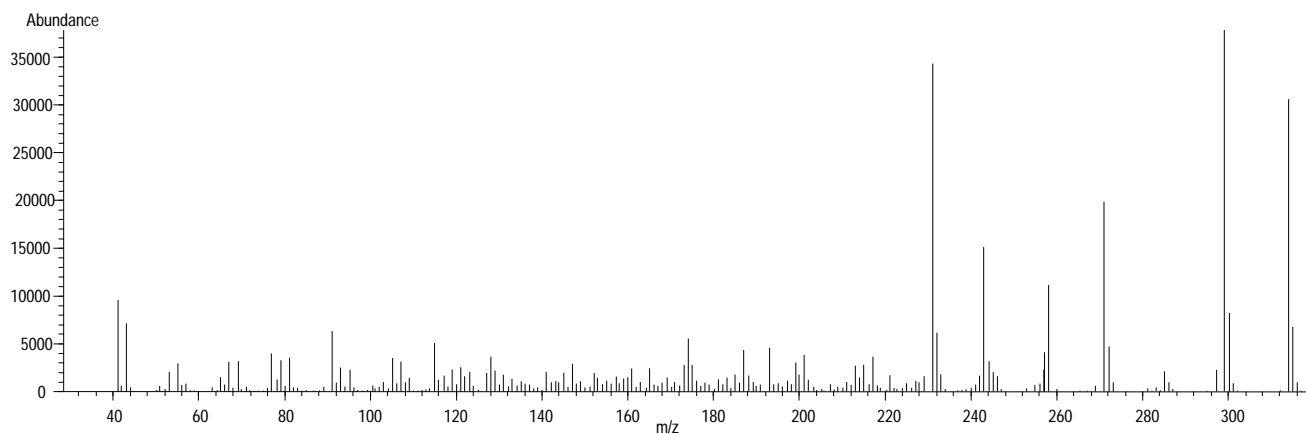
## Capillary Gas Chromatography

**Column** Supelco SPB-1 (Cat. No. 2-4162)  
**I.D.** 0.20 mm **Length** 15 m  
**Film Thickness** 0.20  $\mu$ m  
**Oven Temperature** 50  $^{\circ}$ C (2 min.), then 15 $^{\circ}$ /min. to 300  $^{\circ}$ 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

## 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)

T 4764/93K8811/09/22/03

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.