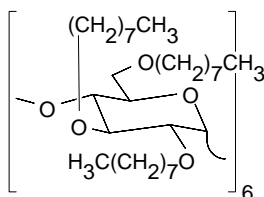

(-)-(1R,2S)-Ephedrine

**Hexakis-(2,3,6-tri-O-octyl)- α -cyclodextrin**(2,3,6-Tri-O-octyl- α -cyclodextrin) $C_{180}H_{348}O_{30}$ M_r 2992.67

[140395-31-9]

[52584](#)**Selectophore[®]**, $\geq 99.0\%$ (TLC)100 mg

Electrochemical Transduction

Ion-Selective Electrodes

Application and Sensor Type ¹

Determination of the enantiomeric purity of ephredrine in the presence of serum cations with solvent polymeric membrane electrode based on Hexakis-(2,3,6-tri-O-octyl)- α -cyclodextrin.

Recommended Membrane Composition

| | | |
|-------|-----|---|
| 2.00 | wt% | Hexakis-(2,3,6-tri-O-octyl)- α -cyclodextrin (52584) |
| 65.60 | wt% | Bis(1-butylpentyl)adipate (02150) |
| 0.40 | wt% | Potassium tetrakis(4-chlorophenyl)borate (60591) |
| 32.00 | wt% | Poly(vinyl chloride) high molecular weight (81392) |

Recommended Cell Assembly

Reference || sample solution || liquid membrane | 0.001 M NH₄Cl | AgCl, Ag

Electrode Characteristics and Function

Selectivity coefficients $\log K_{\text{Eph}^- \text{HCl}, \text{M}}^{\text{Pot}}$ as obtained by the fixed interference method.

$\log K_{(+)-\text{Eph}^- \text{HCl}, (\text{Na}, \text{K}, \text{Ca})}^{\text{Pot}}$ -3.9

$\log K_{(-)-\text{Eph}^- \text{HCl}, (\text{Na}, \text{K}, \text{Ca})}^{\text{Pot}}$ -3.5

¹⁾ Ion background of serum levels

Slope of linear regression: 59 mV at 25°C (ion background of serum levels: 150mM NaCl, 4.3 mM KCl, 1.26 mM CaCl₂)

Enantioselectivity: $\log K_{+/-}^{\text{Pot}}$ -2.7

Detection limit: $\log a_{(+)-\text{Eph}^- \text{HCl}}$ -4.7
 $\log a_{(-)-\text{Eph}^- \text{HCl}}$ -4.4

Lifetime: ~3 months

¹ R. Katakay, P.S. Bates, D. Parker, Functionalized α -cyclodextrins as potentiometric chiral sensors. **Analyst** **117**, 1313 (1992).