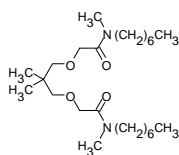


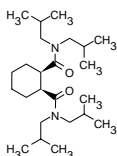
## Lithium



### Lithium ionophore I

(ETH 149; *N,N'*-Diheptyl-*N,N',5,5*-tetramethyl-3,7-dioxanonanediamide)  
 $C_{25}H_{50}N_2O_4$   $M_r = 442.67$  [58821-96-8]

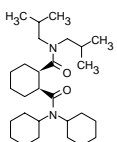
[62557](#) Selectophore®, function tested 50 mg



### Lithium ionophore II

(ETH 1644; *N,N,N',N'*-Tetraisobutyl-*cis*-cyclohexane-1,2-dicarboxamide)  
 $C_{24}H_{46}N_2O_2$   $M_r = 394.64$  [80547-18-8]

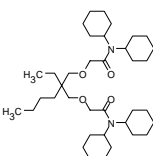
[62559](#) Selectophore®, function tested 50 mg



### Lithium ionophore III

(ETH 1810; *N,N*-Dicyclohexyl-*N',N'*-diiisobutyl-*cis*-cyclohexane-1,2-dicarboxamide)  
 $C_{28}H_{50}N_2O_2$   $M_r = 602.95$  [99281-50-2]

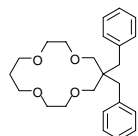
[62558](#) Selectophore®, function tested 50 mg



### Lithium ionophore IV

(ETH 2137; 5-Butyl-5-ethyl-*N,N,N',N'*-tetracyclohexyl-3,7-dioxaazelaic diamide)  
 $C_{37}H_{66}N_2O_4$   $M_r = 602.95$  [108083-23-4]

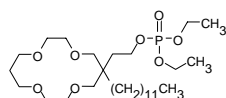
[62561](#) Selectophore®, function tested 50 mg (solution in 0.5 mL heptane)



### Lithium ionophore VI

(6,6-Dibenzyl-14-crown-4; 6,6-Dibenzyl-1,4,8-11-tetraoxacyclotetradecane)  
 $C_{24}H_{32}O_4$   $M_r = 384.52$  [106868-21-7]

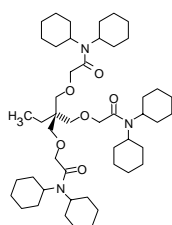
[62567](#) Selectophore®, function tested 10 mg, 100 mg



### Lithium ionophore VII

(6-[2-(Diethylphosphonoxy)ethyl]-6-dodecyl-14-crown-4; 6-Dodecyl(14-crown-4)-6-ethanol diethyl phosphate)  
 $C_{28}H_{57}O_8P$   $M_r = 552.73$  [106868-29-5]

[62569](#) Selectophore®, function tested 10 mg (solution in 0.1 mL heptane)



### Lithium ionophore VIII

(*N,N,N',N',N'',N''*-Hexacyclohexyl-4,4',4''-propylidynetris(3-oxabutryamide))  
 $C_{48}H_{83}N_3O_6$   $M_r = 798.16$  [133338-85-9]

[62571](#) Selectophore®, function tested 50 mg

# Electrochemical Transduction

## Ion-Selective Electrodes

### Application 1 and Sensor Type <sup>1</sup>

Assay of Li<sup>+</sup> activity in biological systems with solvent polymeric membrane electrode based on Lithium ionophore I.

#### Recommended Membrane Composition

1.00	wt%	Lithium ionophore I ( <a href="#">62557</a> )
62.80	wt%	Tris(2-ethylhexyl) phosphate (TEHP) ( <a href="#">93299</a> )
33.00	wt%	Poly(vinyl chloride) high molecular weight ( <a href="#">81392</a> )

#### Recommended Cell Assembly

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

#### Electrode Characteristics and Function

Selectivity Coefficients  $\log K_{Li, M}^{Pot}$  as obtained by the separate solution method (0.1 M solutions of the chlorides)

$\log K_{Li, H}^{Pot}$	-0.1	$\log K_{Li, NH_4}^{Pot}$	-1.3
$\log K_{Li, Na}^{Pot}$	-1.3	$\log K_{Li, Mg}^{Pot}$	-3.7
$\log K_{Li, K}^{Pot}$	-2.1	$\log K_{Li, Ca}^{Pot}$	-3.3

Detection limit:  $\sim 6 \cdot 10^{-6}$  M LiCl

### Application 2 and Sensor Type <sup>2, 3</sup>

Assay of Li<sup>+</sup> activity in serum (therapeutical Li<sup>+</sup> range) with solvent polymeric membrane electrode based on Lithium ionophore II.

#### Recommended Membrane Composition

1.00	wt%	Lithium ionophore II (ETH 1644) ( <a href="#">62559</a> )*
0.40	wt%	Potassium tetrakis( <i>p</i> -chlorophenyl)borate (KTPCIPB) ( <a href="#">60591</a> )
65.60	wt%	2-Nitrophenyl octyl ether ( <a href="#">73732</a> )
33.00	wt%	Poly(vinyl chloride) high molecular weight ( <a href="#">81392</a> )

\* Membranes based on Lithium ionophore III exhibit improved selectivity for measurements in blood.

#### Recommended Cell Assembly

Reference | | sample solution | | liquid membrane | 0.1 M LiCl | AgCl, Ag

#### Electrode Characteristics and Function

Selectivity Coefficients $\log K_{Li, M}^{Pot}$	Required (blood <sup>1)</sup> )	Found
	$\log K_{Li, H}^{Pot}$	<-2.1
$\log K_{Li, Na}^{Pot}$	<-4.3	-2.1
$\log K_{Li, K}^{Pot}$	<-2.8	-2.2
$\log K_{Li, Mg}^{Pot}$	<-3.5	-3.0
$\log K_{Li, Ca}^{Pot}$	<-3.6	-1.7

<sup>1)</sup> for therapeutical Li<sup>+</sup> concentrations (1% interference, worst case)<sup>4, 5</sup>

<sup>1</sup> M. Güggi, U. Fiedler, E. Pretsch, W. Simon, A lithium ion-selective electrode based on a neutral carrier. **Anal. Lett.** **8**, 857 (1975).

<sup>2</sup> D. Erne, D. Ammann, A.F. Zhukov, F. Behm, E. Pretsch, W. Simon, Lipophilic diamides as ionophores for alkali and alkaline earth metal cations. **Helv. Chim. Acta** **65**, 538 (1982).

<sup>3</sup> E. Metzger, D. Ammann, U. Schefer, E. Pretsch, W. Simon, Lipophilic neutral carriers for lithium selective liquid membrane electrodes. **Chimia** **38**, 440 (1984).

**Application 3 and Sensor Type**<sup>3,6,7,8</sup>

 Assay of Li<sup>+</sup> activity in whole blood, plasma and serum (therapeutic Li<sup>+</sup> range) with solvent polymeric membrane electrode based on Lithium ionophore III.

**Recommended Membrane Composition**

1.20	wt%	Lithium ionophore III (ETH 1810) ( <a href="#">62558</a> )
0.40	wt	Potassium tetrakis(4-chlorophenyl)borate ( <a href="#">60591</a> )
65.60	wt%	2-Nitrophenyl octyl ether ( <a href="#">73732</a> )
32.80	wt%	Poly(vinyl chloride) high molecular weight ( <a href="#">81392</a> )

**Recommended Cell Assembly**

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

**Electrode Characteristics and Function**

Selectivity Coefficients $\log K_{Li, M}^{Pot}$	Required (blood <sup>1)</sup> )	Found
$\log K_{Li, H}^{Pot}$	<-2.1	1.0
$\log K_{Li, Na}^{Pot}$	<-4.3	-2.3 (-2.45) <sup>3)</sup>
$\log K_{Li, K}^{Pot}$	<-2.8	-2.6
$\log K_{Li, Mg}^{Pot}$	<-3.5	-4.0
$\log K_{Li, Ca}^{Pot}$	<-3.6	-2.7
Lifetime: $\log P_{TLC}^{(2)}$ ionophore	>8.4	8.3
$\log P_{TLC}^{(2)}$ plasticizer	>12.8	9.3

Stability: Drift [mV/h] 0.02

<sup>1)</sup> for therapeutic Li<sup>+</sup> concentrations (1% interference, worst case)<sup>4,5</sup>
<sup>2)</sup> lipophilicity, determined by thin layer chromatography<sup>9</sup>
<sup>3)</sup> fixed interference method

<sup>4</sup> A. Lewenstam, Ion selective electrodes in clinical chemistry. **Anal. Proc.** **28**, 106 (1991)

<sup>5</sup> U. Oesch, P. Anker, D. Ammann, W. Simon, in: Ion-Selective Electrodes, Eds. E. Pungor, I. Buzás, Akadémiai Kiadó, Budapest 81 (1985).

<sup>6</sup> E. Metzger, D. Ammann, R. Asper, W. Simon, Ion selective liquid membrane electrode for the assay of lithium in blood serum. **Anal. Chem.** **58**, 132 (1986)

<sup>7</sup> E. Metzger, R. Dohner, W. Simon, D.J. Vonderschmitt, K. Gautschi, Lithium/sodium ion concentration ratio measurements in blood serum with lithium and sodium ion selective liquid membrane electrodes. **Anal. Chem.** **59**, 1600 (1987).

<sup>8</sup> J.H. Sim, K.M. Lee, J.S. Lee, D.H. Cho, H. Nam, G.S. Cha, Lithium Ion-Selective Electrode with Improved Lifetime. **Bull. Korean Chem. Soc.** **22**, 765 (2001)

<sup>9</sup> O. Dinten, U.E. Spichiger, N. Chaniotakis, P. Gehrig, B. Rusterholz, W.E. Morf, W. Simon, Lifetime of neutral-carrier-based liquid membranes in aqueous samples and blood and the lipophilicity of membrane components. **Anal. Chem.** **63**, 596 (1991).

#### Application 4 and Sensor Type <sup>7,10</sup>

Determination of Li<sup>+</sup>/Na<sup>+</sup> concentration ratios in blood serum (therapeutic Li<sup>+</sup> range) with solvent polymeric membrane electrodes based on Lithium ionophore IV, combined with a Na<sup>+</sup>-selective polymeric membrane electrode.

##### Recommended Membrane Composition

2.00	wt%	Lithium ionophore IV (ETH 2137) ( <a href="#">62561</a> )
65.60	wt%	Bis(1-butylpentyl) adipate ( <a href="#">02150</a> )
32.40	wt%	Poly(vinyl chloride) high molecular weight ( <a href="#">81392</a> )

##### Recommended composition of the Na<sup>+</sup>-selective solvent polymeric membrane

1.00	wt%	Sodium ionophore III (ETH 2120) ( <a href="#">71734</a> )
66.00	wt%	Bis(1-butylpentyl)decane-1,10-diyl diglutarate (ETH 469) ( <a href="#">30585</a> )
33.00	wt%	Poly(vinyl chloride) high molecular weight ( <a href="#">81392</a> )

##### Recommended Cell Assembly

AgCl,Ag | 0.001 M LiCl, 0.14 M NaCl in 0.5 % aqueous agar | Na<sup>+</sup>-selective membrane | | sample solution | Li<sup>+</sup>-selective membrane liquid membrane | 0.001 M LiCl, 0.14 M NaCl in 0.5% aqueous agar | AgCl,Ag

##### Electrode and System Characteristics

The electrode response function for Li<sup>+</sup>/Na<sup>+</sup> concentration ratio is described by the following equation:

$$E_{ratio} = E_{ratio}^0 + s \cdot \log\left(\frac{C_{Li}}{C_{Na}} + K_{LiNa}^{Pot}\right)$$

Selectivity coefficient  $\log K_{Li,Na}^{Pot} -1.9$

Response time: ~20 s

Residual standard deviation (EMF response function): <0.1 mV

#### Application 5 and Sensor Type <sup>11,12,13</sup>

Neutral carrier for Li<sup>+</sup> in ion-selective PVC membrane electrodes with high selectivity over potassium and sodium ions.

##### Recommended Membrane Composition

1.00	wt%	Lithium ionophore VI ( <a href="#">62567</a> )
0.70	wt%	Potassium tetrakis(4-chlorophenyl)borate ( <a href="#">60591</a> )
70.30	wt%	2-Nitrophenyl octyl ether ( <a href="#">73732</a> )
28.00	wt%	Poly(vinyl chloride) high molecular weight ( <a href="#">81392</a> )

##### Recommended Cell Assembly

Reference | | sample solution | | liquid membrane | 1 M LiCl | AgCl,Ag

##### Electrode Characteristics and Function

Selectivity Coefficients  $\log K_{Li,M}^{Pot}$  determined by the fixed interference method (0.05 M solution for alkali metal ions and H<sup>+</sup>, 0.5 M for alkaline earth metal ions and NH<sub>4</sub><sup>+</sup>).

$\log K_{Li,H}^{Pot}$	-3.0	$\log K_{Li,NH_4}^{Pot}$	-3.0
$\log K_{Li,Na}^{Pot}$	-2.4	$\log K_{Li,Mg}^{Pot}$	-4.3
$\log K_{Li,K}^{Pot}$	-2.3	$\log K_{Li,Ca}^{Pot}$	-4.7

Slope of linear regression: 59-60 mV/decade

<sup>10</sup> E. Metzger, R. Aeschmann, M. Egli, G. Suter, R. Dohner, D. Ammann, M. Dobler, W. Simon, 3,7-Dioxaazelaamides as Ionophores for Lithium Ion Selective Liquid Membrane Electrodes.. *Helv. Chim. Acta* **69**, 1821 (1986).

<sup>11</sup> K. Kimura, H. Yano, S. Kitazawa, T. Shono, Synthesis and Selectivity for Lithium of Lipophilic 14-Crown-4 Derivatives Bearing Bulky Substituents or an Additional Binding Site in the Side Arm, *J. Chem. Soc. Perkin Trans. II*, 1945 (1986).

<sup>12</sup> K. Kimura, O. Oishi, T. Miura, T. Shono, Lithium Ion Selective Electrodes Based on Crown Ethers for Serum Lithium Assay. *Anal. Chem.* **59**, 2331 (1987).

<sup>13</sup> S. Sawada, T. Osakai, M. Senda, Voltammetric Lithium Ion-Selective Electrodes Based on Ion Transfer at the Oil/Water Interface Facilitated by Neutral Ionophores. *Anal. Sci.* **11**, 733 (1995).

### Application 6 and Sensor Type<sup>14</sup>

Neutral lithium ionophore in ion-selective polymer membrane electrodes with excellent selectivity over potassium and sodium ions.

#### Recommended Membrane Composition

1.00	wt%	Lithium ionophore VII ( <a href="#">62569</a> )
0.70	wt%	Potassium tetrakis(4-chlorophenyl)borate ( <a href="#">60591</a> )
70.20	wt%	2-Nitrophenyl octyl ether ( <a href="#">73732</a> )
28.10	wt%	Poly(vinyl chloride) high molecular weight ( <a href="#">81392</a> )

#### Recommended Cell Assembly

Reference | | sample solution | | liquid membrane | 1 M LiCl | AgCl, Ag

#### Electrode Characteristics and Function

Selectivity Coefficients  $\log K_{Li, M}^{Pot}$  determined by the fixed interference method (0.05 M solution for alkali metal ions and H<sup>+</sup>, 0.5 M for alkaline earth metal ions and NH<sub>4</sub><sup>+</sup>).

$\log K_{Li, H}^{Pot}$	-3.2	$\log K_{Li, NH_4}^{Pot}$	-3.0
$\log K_{Li, Na}^{Pot}$	-2.3	$\log K_{Li, Mg}^{Pot}$	-4.4
$\log K_{Li, K}^{Pot}$	-2.5	$\log K_{Li, Ca}^{Pot}$	-4.0

Slope of linear regression: 59-60 mV

### Application 7 and Sensor Type<sup>15</sup>

Sensor with increased stability and lifetime for assay of Li<sup>+</sup> based on Lithium ionophore VIII.

#### Recommended Membrane Composition

2.50	wt%	Lithium ionophore VIII ( <a href="#">62571</a> )
0.23	wt%	Potassium tetrakis(4-chlorophenyl)borate ( <a href="#">60591</a> )
64.50	wt%	Bis(1-butylpentyl) adipate ( <a href="#">02150</a> )
32.80	wt%	Poly(vinyl chloride) high molecular weight ( <a href="#">81392</a> )

#### Recommended Cell Assembly

Reference | | sample solution | | liquid membrane | 0.01 M LiCl | AgCl, Ag

#### Electrode Characteristics and Function

Selectivity Coefficients  $\log K_{Li, M}^{Pot}$  determined by the separate solution method (0.1 M solutions of the chlorides).

$\log K_{Li, Na}^{Pot}$	-2.4	$\log K_{Li, Mg}^{Pot}$	-4.3
$\log K_{Li, K}^{Pot}$	-2.3	$\log K_{Li, Ca}^{Pot}$	-4.7
$\log K_{Li, NH_4}^{Pot}$	-3.0		

Slope of linear regression: 59-60 mV

<sup>14</sup> K. Kimura, H. Yano, S. Kitazawa, T. Shono, Synthesis and selectivity for lithium of lipophilic 14-crown-4 derivatives bearing bulky substituents or an additional binding site in the side arm. **J. Chem. Soc. Perkin Trans. 2**, 1945 (1986).

<sup>15</sup> M. Bochenska, W. Simon, Performance of a highly lipophilic Li<sup>+</sup> ionophore in solvent polymeric membranes. **Microchim. Acta III**, 277 (1990).