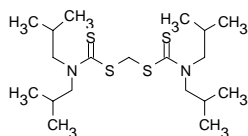


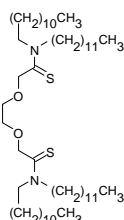
Lead



Lead ionophore II

(MBDiBDTC; S,S'-Methylenebis(N,N-diisobutylthioacetamide)
 $C_{19}H_{38}N_2S_4$ $M_r = 422.77$ [90276-58-7]

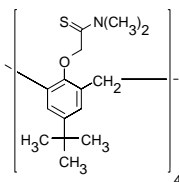
[15336](#) Selectophore®, function tested 50 mg



Lead ionophore III

(ETH 5435; N,N,N',N'-Tetradodecyl-3,6-dioxaocanedithioamide)
 $C_{54}H_{108}N_2O_2S_4$ $M_r = 881.35$ [141754-61-2]

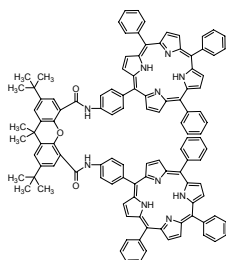
[98108](#) Selectophore®, function tested 50 mg



Lead ionophore IV

(*tert*-Butylcalix[4]arene-tetrakis(N,N-dimethylthioacetamide))
 $C_{60}H_{84}N_4O_4S_4$ $M_r = 1054.59$ [145237-46-3]

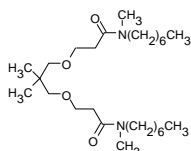
[15343](#) Selectophore®, function tested 50 mg



Lead ionophore VII

(ADPX)
 $C_{113}H_{88}N_{10}O_3$ $M_r = 1633.97$

[61796](#) Selectophore®, function tested 50 mg



Uranyl ionophore I

(ETH 295; N,N'-Diheptyl-N,N',6,6-tetramethyl-4,8-dioxaundecanediamide)
 $C_{27}H_{54}N_2O_4$ $M_r = 470.73$ [69844-41-3]

[94265](#) Selectophore®, function tested 50 mg

Electrochemical Transduction

- Ion-Selective Electrodes
- Chemically modified Electrodes

Optical Transduction

Electrochemical Transduction

Ion-Selective Electrodes

Application 1 and Sensor Type ^{1,2}

Lead-selective solvent polymeric membrane electrode based on Lead ionophore II.

Recommended Membrane Composition

11.20	wt%	Lead ionophore II (15336)
2.00	wt%	Potassium tetrakis(<i>p</i> -chlorophenyl)borate (KTpCIPB) (60591)
49.60	wt%	2-Nitrophenyl octyl ether (73732)
37.20	wt%	Poly(vinyl chloride) high molecular weight (81392)

Recommended Cell Assembly

Reference | | sample solution | | liquid membrane | 0.001 M PbCl₂ | AgCl, Ag

Electrode Characteristics and Function

Selectivity coefficients $\log K_{Pb, M}^{Pot}$ as obtained by the mixed solution method (concentration of the interfering ions are 0.1 M except for Cu (10⁻⁵ M), Fe and Zn (10⁻³ M). Measured in Pb(NO₃)₂ solution of 10⁻¹ to 10⁻⁸ M).

$\log K_{Pb, Na}^{Pot}$	-2.14	$\log K_{Pb, Mn}^{Pot}$	-5.21
$\log K_{Pb, K}^{Pot}$	-2.20	$\log K_{Pb, Ca}^{Pot}$	-5.45
$\log K_{Pb, Fe}^{Pot}$	-2.54	$\log K_{Pb, Mg}^{Pot}$	-5.26
$\log K_{Pb, Zn}^{Pot}$	-3.50	$\log K_{Pb, Sr}^{Pot}$	-5.25
$\log K_{Pb, Cd}^{Pot}$	-3.57	$\log K_{Pb, Cr}^{Pot}$	-4.6
$\log K_{Pb, Ni}^{Pot}$	-5.0	$\log K_{Pb, Cu}^{Pot}$	0.7
$\log K_{Pb, Co}^{Pot}$	-5.2		

Slope of linear regression: 28 mV (10⁻⁶ – 10⁻² M Pb(NO₃)₂).

Detection limit: 3.5·10⁻⁷ M

Response time: 95% response time 16s (10⁻³ to 10⁻² M Pb(NO₃)₂).

pH range: 3.1-5.4 (measured in 10⁻³ M Pb(NO₃)₂ solution).

Application 2 and Sensor Type ³

Lead-selective solvent polymeric membrane electrode based on Lead ionophore III.

Recommended Membrane Composition

1.31	wt%	Lead ionophore III (ETH 5435) (98108)
0.82	wt%	Potassium tetrakis[3,5-bis(trifluoromethyl)phenyl]borate (KTpCIPB) (60588)
62.57	wt%	Bis(2-ethylhexyl) sebacate (84818)
35.30	wt%	Poly(vinyl chloride) high molecular weight (81392)

Recommended Cell Assembly

Reference | | sample solution | | liquid membrane | 0.0005 M PbCl₂ and 0.05M MgCl₂ | AgCl, Ag

¹ S. Kamata, K. Onoyama, Lead-selective membrane electrode using methylene bis(diisobutyldithiocarbamate) neutral carrier. **Anal. Chem.** **63**, 1295 (1991).

² S. Kamata, K. Onoyama, Methylene Bis(diisobutyldithiocarbamate) Neutral Carrier as Lead Sensing Material. **Chem. Lett.** **653** (1991).

³ T. Sokalski, A. Ceresa, T. Zwicky, E. Pretsch, Large Improvement of Lower Detection Limit of Ion-Selective Polymer Membrane Electrodes. **J. Am. Chem. Soc.** **119**, 11347 (1997).

Electrode Characteristics and Function

Selectivity coefficients $\log K_{Pb, M}^{Pot}$ as obtained by the separate solution method in 0.01M nitrate solutions of Na^+ , K^+ , Ca^{2+} , Mg^{2+} , Mn^{2+} , Co^{2+} , Ni^{2+} , Mn^{2+} , Cu^{2+} , Zn^{2+} and Cd^{2+} buffered to pH 4.7 with a magnesium acetate buffer ($2.5 \cdot 10^{-4}$ M Mg^{2+} , 10^{-3} M total Ac)

$\log K_{Pb, H}^{Pot}$	-2.2	$\log K_{Pb, Sr}^{Pot}$	-4.2
$\log K_{Pb, Li}^{Pot}$	-2.7	$\log K_{Pb, Ba}^{Pot}$	-4.1
$\log K_{Pb, Na}^{Pot}$	-2.2	$\log K_{Pb, Mn}^{Pot}$	-4.1
$\log K_{Pb, K}^{Pot}$	-2.2	$\log K_{Pb, Co}^{Pot}$	-4.3
$\log K_{Pb, NH_4}^{Pot}$	-2.5	$\log K_{Pb, Ni}^{Pot}$	-4.2
$\log K_{Pb, Mg}^{Pot}$	-3.6	$\log K_{Pb, Cu}^{Pot}$	-0.5
$\log K_{Pb, Ca}^{Pot}$	-4.0	$\log K_{Pb, Zn}^{Pot}$	-3.7
$\log K_{Pb, Cd}^{Pot}$	0.9		

Application 3 and Sensor Type ^{4, 5}

Lead-selective solvent polymeric membrane electrode based on Lead ionophore V with very good selectivity over Cu, Zn and Cd.

Recommended Membrane Composition

1.00	wt%	Lead ionophore V (15343)
0.35	wt%	Potassium tetrakis(4-chlorophenyl)borate (60591)
65.65	wt%	2-Nitrophenyl octyl ether (73732)
33.00	wt%	Poly(vinyl chloride) high molecular weight (81392)

Recommended Cell Assembly

Reference (e.g. Ag, AgCl, 0.1 M KCl | 0.1M KNO₃) || sample solution || liquid membrane | 0.005 M PbCl₂ | AgCl, Ag

Electrode Characteristics and Function

Selectivity coefficients $\log K_{Pb, M}^{Pot}$ as obtained by the fixed interference method. For Ag and Hg the separate solution method is used (0.01 M solution, pH 4 for Pb and Ag, pH 2 for Hg)

$\log K_{Pb, Hg}^{Pot}$	0.6	$\log K_{Pb, K}^{Pot}$	-4.3
$\log K_{Pb, Ag}^{Pot}$	1.5	$\log K_{Pb, Ba}^{Pot}$	-4.8
$\log K_{Pb, Cu}^{Pot}$	-3.3	$\log K_{Pb, Ca}^{Pot}$	-4.8
$\log K_{Pb, Na}^{Pot}$	-3.5	$\log K_{Pb, Zn}^{Pot}$	-4.8
$\log K_{Pb, Li}^{Pot}$	-3.5	$\log K_{Pb, Mg}^{Pot}$	-5.0

⁴ E. Malinowska, Z. Brzozka, K. Kasiura, R.J.M. Egberink, D.N. Reinhoudt, Lead selective electrodes based on thioamide functionalized calix[4]arenes as ionophores. **Anal. Chim. Acta** **298**, 253 (1994).

⁵ C. McGraw, T. Radu, A. Radu, D. Diamond, Evaluation of Liquid- and Solid-Contact, Pb²⁺-Selective Polymer-Membrane Electrodes for Soil Analysis, **Electroanalysis** **20**, 340 (2008).

$\log K_{\text{Pb, Cd}}^{\text{Pot}}$	-3.7	$\log K_{\text{Pb, Ni}}^{\text{Pot}}$	-5.0
$\log K_{\text{Pb, NH}_4}^{\text{Pot}}$	-4.0	$\log K_{\text{Pb, Co}}^{\text{Pot}}$	-5.0

Slope of linear regression: 28.7 mV/dec ($1 \cdot 10^{-6} - 2 \cdot 10^{-2}$ M Pb^{2+})

Response time: 95% response time: <10 s

Detection limit: $3 \cdot 10^{-7}$ Pb^{2+}

pH range: 3-6

Drift (0.01 M Pb, t = 25 min): <0.01 mV/min

Application 4 and Sensor Type ⁶

Lead-selective solvent polymeric membrane electrode based on Lead ionophore VII at pH 4.5-7.5.

Recommended Membrane Composition

3.0	wt%	Lead ionophore VII (61796)
2.8	wt%	Sodium tetrphenylborate (NaTPB) (72018)
63.2	wt%	2-Nitrophenyl octyl ether (o-NPOE) (73732)
31.0	wt%	Poly(vinyl chloride) high molecular weight (81392)

Recommended Cell Assembly

Reference (e.g. Ag, AgCl, KCl (sat) | sat. NH_4NO_3) || sample solution || liquid membrane | 0.001 M $\text{Pb}(\text{NO}_3)_2$ | AgCl, Ag

Electrode Characteristics and Function

Selectivity coefficients $\log K_{\text{Pb, M}}^{\text{Pot}}$ as obtained by the separate solution method. (0.01M solutions in 0.001 M Tris- HNO_3 buffer, pH 6.0)

$\log K_{\text{Pb, Na}}^{\text{Pot}}$	-1.0	$\log K_{\text{Pb, Mg}}^{\text{Pot}}$	-3.3
$\log K_{\text{Pb, K}}^{\text{Pot}}$	-0.6	$\log K_{\text{Pb, Zn}}^{\text{Pot}}$	-2.8
$\log K_{\text{Pb, Ca}}^{\text{Pot}}$	-3.2	$\log K_{\text{Pb, Cu}}^{\text{Pot}}$	-14.5

Slope of linear regression: 28.4 mV/dec ($4 \cdot 10^{-5} - 5 \cdot 10^{-2}$ M Pb^{2+})

Detection limit: $1 \cdot 10^{-5}$ Pb^{2+}

Application 5 and Sensor Type ⁷

Lead-selective solvent polymeric membrane electrode based on Uranyl ionophore I.

Recommended Membrane Composition

1.00	wt%	Uranyl ionophore I (94265)
0.42	wt%	Potassium tetrakis(4-chlorophenyl)borate (60591)
68.58	wt%	2-Nitrophenyl octyl ether (73732)
30.00	wt%	Poly(vinyl chloride) high molecular weight (81392)

Recommended Membrane Composition

Reference (e.g. Ag, AgCl, 1 M KCl | 0.1 M KNO_3) | sample solution | liquid membrane | 0.005 M PbCl_2 | AgCl, Ag

⁶ W.J. Zhang, C.Y. Li, Y.B. Zhang, J. Jin, Synthesis of an Amide-Linked Diporphyrin Xanthene as a Neutral Carrier for a Lead(II)-Sensitive Electrode. **Anal. Lett.** **40**, 1023 (2007).

⁷ E. Malinowska, Lead-selective membrane electrodes based on neutral carriers. Part I. Acyclic amides and oxamides. **Analyst** **115**, 1085 (1990).

Electrode Characteristics and Function

Selectivity coefficients $\log K_{Pb, M}^{Pot}$ as obtained by the separate solution method (pH 4).

$\log K_{Pb, Ag}^{Pot}$	0.16	$\log K_{Pb, Ca}^{Pot}$	-2.3
$\log K_{Pb, H}^{Pot}$	-0.16	$\log K_{Pb, Cu}^{Pot}$	-3.9
$\log K_{Pb, Na}^{Pot}$	-3.6	$\log K_{Pb, Cd}^{Pot}$	-4.1
$\log K_{Pb, Li}^{Pot}$	-2.9	$\log K_{Pb, Zn}^{Pot}$	-4.3
$\log K_{Pb, NH_4}^{Pot}$	-3.5	$\log K_{Pb, Mg}^{Pot}$	-4.6
$\log K_{Pb, K}^{Pot}$	-3.9	$\log K_{Pb, Co}^{Pot}$	-4.1
$\log K_{Pb, Sr}^{Pot}$	-1.7	$\log K_{Pb, Ni}^{Pot}$	-4.6
$\log K_{Pb, Ba}^{Pot}$	-2.4		

Slope of linear regression: 35.3 mV ($6 \cdot 10^{-6}$ - $1 \cdot 10^{-1}$ M Pb^{2+})

Detection limit: $3 \cdot 10^{-6}$ M $Pb(NO_3)_2$

Response time: 95% response time: <15 s (10^{-3} to 10^{-2} M Pb^{2+})

pH range: 2-6

Drift: <0.01 mV/min (10^{-2} M Pb^{2+})

Chemically modified Electrodes

Application and Sensor Type⁸

Lead-selective glassy carbon electrodes for accumulation voltametry based on Lead ionophore II. With this chemically modified electrode (CME) mmol concentration of lead(II) can be detected.

Recommended Membrane Composition

Solution A: $5 \cdot 10^{-3}$ M Lead ionophore II ([15336](#)) in 2-propanol ([34863](#))

Solution B: Nafion (5% solution in lower aliphatic alcohols and water) ([70160](#)) diluted to 1% with methanol ([42105](#))

Coating solutions:

2 μ L Solution A

2 μ L Solution B

10 μ L Methanol ([42105](#))

A glassy carbon electrode is polished on Al_2O_3 slurry, sonicated in methanol, air-dried and coated with the solution at an area of 0.38 cm^2 . This corresponds to a coating with 10 mmol Lead ionophore II and 20 μ g Nafion with a calculated film thickness of 0.39 μ m.

Electrode Characteristics and Function

Accumulation of lead(II) from 0.1 M acetate buffer (pH 4) followed by voltametry gave a quasi-reversible reduction peak at -0.85 V.

Half-wave potential: $E_{1/2}$: -0.57 V

Peak to peak separation: $\Delta E_p = 0.51$ V

Linear response range: $5 \cdot 10^{-7}$ M to $6 \cdot 10^{-6}$ M Pb^{2+} (120 s accumulation)

Detection limit: 10^{-7} M Pb^{2+} (with 600s accumulation)

Interference: Ca^{2+} , Mg^{2+} , Ni^{2+} , Cu^{2+} , Cd^{2+} , Ag^+ and Hg^{2+} are the most highly interfering species.

⁸ M. Lerchi, E. Bakker, B. Rusterholz, W. Simon, Lead-selective bulk optodes based on neutral ionophores with subnanomolar detection limits. *Anal. Chem.* **64**,1534 (1992).

Optical Transduction

Application and Sensor Type^{9,10}

Lead-selective bulk optode membrane with subnanomolar detection limit based on Lead ionophore III.

Recommended Membrane Composition

2.40	wt%	Lead ionophore III (98108)
0.72	wt%	Chromoionophore VII (ETH 5418) (27095)
64.00	wt%	Bis(2-ethylhexyl) sebacate (84818)
0.88	wt%	Potassium tetrakis[3,5-bis(trifluoromethyl)phenyl] borate (60588)
32.00	wt%	Poly(vinyl chloride) high molecular weight (81392)

Optode Characteristics and Function

Selectivity coefficients $\log K_{\text{Pb, M}}^{\text{Opt}}$ as obtained by the separate solution method (pH 4.7).

$\log K_{\text{Pb, Na}}^{\text{Opt}}$	-5.6	$\log K_{\text{Pb, Co}}^{\text{Opt}}$	-6.5
$\log K_{\text{Pb, K}}^{\text{Opt}}$	-5.3	$\log K_{\text{Pb, Ni}}^{\text{Opt}}$	-4.8
$\log K_{\text{Pb, Mg}}^{\text{Opt}}$	-10.9	$\log K_{\text{Pb, Cu}}^{\text{Opt}}$	0.7
$\log K_{\text{Pb, Ca}}^{\text{Opt}}$	-10.8	$\log K_{\text{Pb, Zn}}^{\text{Opt}}$	-4.6
$\log K_{\text{Pb, Mn}}^{\text{Opt}}$	-3.0	$\log K_{\text{Pb, Cd}}^{\text{Opt}}$	1.2

Measuring range: $5 \cdot 10^{-9}$ to $5 \cdot 10^{-3}$ M Pb^{2+} (pH 5.68)

Detection limit: $3.2 \cdot 10^{-12}$ M Pb^{2+}

⁹ M. Lerchi, W. Simon, Optodes for environmental chemical monitoring of lead. *Proc. SPIE Int. Soc. Opt. Eng.* **1716**, 336 (1993).

¹⁰ E. Bakker, M. Willer, E. Pretsch, Detection limit of ion-selective bulk optodes and corresponding electrodes. *Anal. Chim. Acta* **282**, 265 (1993).