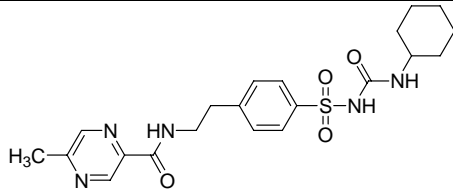


Samarium

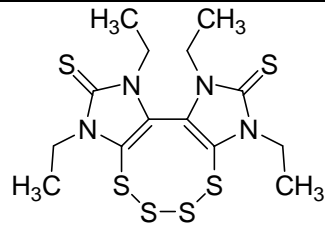


Samarium ionophore I

(Glipizide; 1-Cyclohexyl-3-{4-[2-(5-methylpyrazine-2-carboxamido)ethyl]phenylsulfonyl}urea)

$C_{21}H_{27}N_5O_4S$ M_r 445.54 [29094-61-9]

[30553](#) **Selectophore®** 25 mg, 100 mg



Samarium ionophore II

(1,3,8,10-Tetraethyl-1,3,8,10-tetrahydro[1,2,3,4]tetrathioino[5,6-d:7,8-d']diimidazole-2,9-dithione; Et₄todit)

$C_{14}H_{20}N_4S_6$ M_r 436.73 [120097-53-2]

[15906](#) **Selectophore®** 50 mg

Electrochemical Transduction

Ion-Selective Electrodes

Application 1 and Sensor Type ¹

Assay of Sm^{3+} activity in aqueous solution with solvent polymeric membrane electrode based on Samarium ionophore I.

Recommended Membrane Composition

11	wt%	Samarium ionophore I (30553)
55	wt%	Benzyl acetate (43957)
4	wt%	Sodium tetraphenylborate (NaTPB) (72018)
30	wt%	Poly(vinyl chloride) high molecular weight (81392)

Recommended Cell Assembly

Reference || sample solution || liquid membrane | 0.001 M SmCl_3 | AgCl, Ag

Electrode Characteristics and Function

Selectivity coefficients $\log K_{\text{Sm, M}}^{\text{Pot}}$ as obtained by the matched potential method. Optimum conditioning time for the membrane sensor in a $1.0 \cdot 10^{-3}$ M SmCl_3 solution is 12 h.

$\log K_{\text{Sm, La}}^{\text{Pot}}$	-2.4	$\log K_{\text{Sm, Co}}^{\text{Pot}}$	-3.1
$\log K_{\text{Sm, Ce}}^{\text{Pot}}$	-2.1	$\log K_{\text{Sm, Ni}}^{\text{Pot}}$	-2.7
$\log K_{\text{Sm, Gd}}^{\text{Pot}}$	-2.4	$\log K_{\text{Sm, Sr}}^{\text{Pot}}$	-2.6
$\log K_{\text{Sm, Pb}}^{\text{Pot}}$	-2.0	$\log K_{\text{Sm, Ca}}^{\text{Pot}}$	-2.7
$\log K_{\text{Sm, Hg}}^{\text{Pot}}$	-2.7	$\log K_{\text{Sm, Mg}}^{\text{Pot}}$	-2.6
$\log K_{\text{Sm, Ag}}^{\text{Pot}}$	-2.0	$\log K_{\text{Sm, Na}}^{\text{Pot}}$	-3.1
$\log K_{\text{Sm, Cu}}^{\text{Pot}}$	-2.8	$\log K_{\text{Sm, K}}^{\text{Pot}}$	-3.1

Slope of linear regression: 19.8 mV/dec ($1.0 \cdot 10^{-6}$ to $1.0 \cdot 10^{-1}$ M SmCl_3)

Detection limit: $6.0 \cdot 10^{-7}$ M Sm^{3+}

Response time: 95% response time: <10 s ($>10^{-5}$ M Sm^{3+})

Lifetime: >2 months

¹ M.R. Ganjali, M.R. Pourjavid, M. Rezapour, S. Haghgoo, Novel Samarium(III) selective membrane sensor based on glipizid. **Sensors and Actuators B 89**, 21 (2003).

Application 2 and Sensor Type ²

Assay of Sm³⁺ activity in aqueous solution with solvent polymeric membrane electrode based on Samarium ionophore II.

Recommended Membrane Composition

3	wt%	Samarium ionophore II (15906)
2	wt%	Sodium tetraphenylborate (NaTPB) (72018)
65	wt%	2-Nitrophenyl octyl ether (73732)
30	wt%	Poly(vinyl chloride) high molecular weight (81392)

Recommended Cell Assembly

Reference	sample solution	liquid membrane	0.001 M Sm(NO ₃) ₃	KCl (3M)	AgCl, Ag	(PME)
Reference	sample solution	liquid membrane	graphite surface			(CGE)

Electrode Characteristics and Function

Selectivity coefficients $\log K_{Sm, M}^{Pot}$ as obtained by the matched potential method. Optimum conditioning time for the membrane sensor in a $1.0 \cdot 10^{-3}$ M Sm(NO₃)₃ solution is 24 h.²

	PME	CGE		PME	CGE
$\log K_{Sm, La}^{Pot}$	-1.3	-3.2	$\log K_{Sm, Co}^{Pot}$	-2.3	-3.3
$\log K_{Sm, Ce}^{Pot}$	-1.1	-3.1	$\log K_{Sm, Ni}^{Pot}$	-2.4	-3.4
$\log K_{Sm, Gd}^{Pot}$	-1.1	-3.1	$\log K_{Sm, Tl}^{Pot}$	-3.2	-4.3
$\log K_{Sm, Pb}^{Pot}$	-1.0	-2.2	$\log K_{Sm, Ca}^{Pot}$	-2.9	-3.4
$\log K_{Sm, Hg}^{Pot}$	-1.1	-2.2	$\log K_{Sm, Mg}^{Pot}$	-2.7	-3.6
$\log K_{Sm, Ag}^{Pot}$	-2.4	-3.4	$\log K_{Sm, Ba}^{Pot}$	-2.7	-3.5
$\log K_{Sm, Cu}^{Pot}$	-2.5	-3.4	$\log K_{Sm, Na}^{Pot}$	-3.5	-4.1
$\log K_{Sm, Cd}^{Pot}$	-2.7	-3.7	$\log K_{Sm, K}^{Pot}$	-3.4	-4.6
$\log K_{Sm, Zn}^{Pot}$	-2.1	-3.9	$\log K_{Sm, Li}^{Pot}$	-3.4	-4.8

PME (Polymeric Membrane Electrode):

Conditions	Reference solution: $1.0 \cdot 10^{-6}$ M Sm ³⁺ : $1.0 \cdot 10^{-2}$ M - $1.0 \cdot 10^{-5}$ M M ⁿ⁺ : $1.0 \cdot 10^{-2}$ M - $1.0 \cdot 10^{-3}$ M
------------	--

Slope of linear regression:	19.6 ± 0.5 mV/dec ($1.0 \cdot 10^{-5}$ to $1.0 \cdot 10^{-1}$ M Sm(NO ₃) ₃)
Detection limit:	$8.0 \cdot 10^{-6}$ M Sm ³⁺
Response time:	20 s
pH range:	4.0-6.5

CGE (Coated Graphite Electrode):

Conditions:	Reference solution: $1.0 \cdot 10^{-8}$ M Sm ³⁺ : $1.0 \cdot 10^{-3}$ M - $1.0 \cdot 10^{-7}$ M M ⁿ⁺ : $1.0 \cdot 10^{-2}$ M - $1.0 \cdot 10^{-4}$ M
Slope of linear regression:	19.3 ± 0.3 mV/dec ($1.0 \cdot 10^{-7}$ to $1.0 \cdot 10^{-1}$ M Sm(NO ₃) ₃)
Detection limit:	$1.6 \cdot 10^{-9}$ M Sm ³⁺
Response time:	18 s

² M. Shamsipur, M. Hosseini, K. Alisadeh, Z. Talebpour, M.F. Mousavi, M.R. Ganjali, M. Arca, V. Lippolis, PVC Membrane and Coated Graphite Potentiometric Sensors Based on Et₄todit for Selective Determination of Samarium(III). **Anal. Chem** **75**, 5680 (2003).