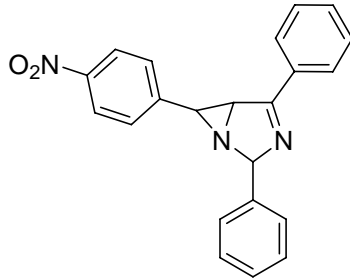


Tin



Tin ionophore II

(6-(4-Nitrophenyl)-2,4-diphenyl-1,3-diazabicyclo[3.1.0]hex-3-ene)
C₂₂H₁₇N₃O₂ M_r 355.39 [13591-58-7]

[73443](#) **Selectophore[®], function tested** 50 mg

Electrochemical Transduction

Ion-Selective Electrodes

Application 1 and Sensor Type ¹

Assay of Sn²⁺ activity in aqueous solution with solvent polymeric membrane electrode based on Tin ionophore II.

Recommended Membrane Composition

5.5	wt%	Tin ionophore II (73443)
58.0	wt%	Acetophenone (00790)
6.5	wt%	Oleic acid (05508)
30.0	wt%	Poly(vinyl chloride) high molecular weight (81392)

Recommended Cell Assembly

Reference || sample solution || liquid membrane | 0.001 M SnCl₂ + 0.001 M HCl | AgCl, Ag

Electrode Characteristics and Function

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

$\log K_{Sn, Ca}^{Pot}$	-1.0	$\log K_{Sn, Zn}^{Pot}$	-1.0
$\log K_{Sn, Mg}^{Pot}$	-1.3	$\log K_{Sn, Ni}^{Pot}$	-3.1
$\log K_{Sn, Co}^{Pot}$	-1.0	$\log K_{Sn, Cu}^{Pot}$	-1.3

Slope of linear regression: 30.1 mV ($4 \cdot 10^{-10}$ to 0.033 mol/L)

Detection limit: $2 \cdot 10^{-10}$ M Zn²⁺

¹ M. Arvand, A.M. Moghimi, A. Afshari, N. Mahmoodi, Potentiometric membrane sensor based on 6-(4-nitrophenyl)-2,4-diphenyl-3,5-diaza-bicyclo[3.1.0]hex-2-ene for detection of Sn(II) in real samples. **Anal. Chim. Acta** **579**, 102 (2006).