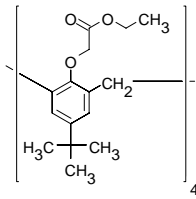


Aldehyde



Sodium ionophore X

(4-*tert*-Butylcalix[4]arene-tetraacetic acid tetraethyl ester)
 $C_{60}H_{80}O_{12}$ M_r 993.29 [97600-39-0]

[71747](#) **Selectophore®** 50 mg

Electrochemical Transduction

Ion-Selective Electrodes

Application 1 and Sensor Type ¹

Aldehyde(heptanal)-selective solvent polymeric membrane electrode based on Sodium ionophore X. The aldehyde is reacted in situ with [Girard's reagent P](#) to generate the hydrazone which is recognised by the electrode.

Recommended Membrane Composition

4.50	wt%	Sodium ionophore X (71747)
60.30	wt%	Bis(1-ethylhexyl) phthalate (80030)
35.20	wt%	Poly(vinyl chloride) high molecular weight (81392)

Recommended Cell Assembly

Reference || sample solution || liquid membrane | 10⁻⁴ M heptanal-Girard's reagent adduct | AgCl, Ag

Electrode Characteristics and Function

Selectivity coefficients $\log K_{\text{Heptanal, X}}^{\text{Pot}}$ as obtained by the separate solution method (solution with 0.01 M ammonium acetate as supporting electrolyte (pH 6). Concentration of interfering species (except NH₄) 2•10⁻⁴ mol/L).

$\log K_{\text{Heptanal, HCHO}}^{\text{Pot}}$	-1.8	$\log K_{\text{Heptanal, NH}_4}^{\text{Pot}}$	-5.4
$\log K_{\text{Heptanal, Butanal}}^{\text{Pot}}$	-1.7	$\log K_{\text{Heptanal, K}}^{\text{Pot}}$	-2.8
$\log K_{\text{Heptanal, Vanilin}}^{\text{Pot}}$	-1.6	$\log K_{\text{Heptanal, Na}}^{\text{Pot}}$	-0.3
$\log K_{\text{Heptanal, n - Butylamine}}^{\text{Pot}}$	-1.9		

Slope of linear regression: 68 ± 1 mV (10⁻⁶ – 10⁻² M heptanal)

Detection limit: 9•10⁻⁷ mol/L heptanal

Application 2 and Sensor Type ²

Formaldehyde-selective solvent polymeric membrane electrode based on Sodium ionophore X. The aldehyde is reacted in situ with a modified Girard's reagent G2 to generate a lipophilic hydrazone which is recognised by the electrode.

Recommended Membrane Composition

4.50	wt%	Sodium ionophore X (71747)
60.30	wt%	Bis(1-ethylhexyl) phthalate (80030)
35.20	wt%	Poly(vinyl chloride) high molecular weight (81392)

Recommended Cell Assembly

Reference || sample solution || liquid membrane | 10⁻³ M heptanal (modified)-Girard's reagent adduct | AgCl, Ag

Electrode Characteristics and Function

Selectivity coefficients $\log K_{\text{Heptanal, X}}^{\text{Pot}}$ as obtained by the separate solution method (solution with 0.01 M ammonium acetate as supporting electrolyte (pH 5.4)).

$\log K_{\text{Heptanal, n - Propanol}}^{\text{Pot}}$	-3.6	$\log K_{\text{Heptanal, NH}_4}^{\text{Pot}}$	-3.5
$\log K_{\text{Heptanal, n - Butylamine}}^{\text{Pot}}$	-1.1	$\log K_{\text{Heptanal, K}}^{\text{Pot}}$	-2.6
$\log K_{\text{Heptanal, Butanal}}^{\text{Pot}}$	0.8	$\log K_{\text{Heptanal, Na}}^{\text{Pot}}$	0.9

Slope of linear regression: 32.4 mV (4•10⁻⁵ – 10⁻² M formaldehyde)

Detection limit: 1.2•10⁻⁵ mol/L formaldehyde.

¹ W.H. Chan, P. X. Cai, X. H. Gu, Aldehyde-selective polymeric membrane electrodes based on a calix[4]arene ionophore. **Analyst** **119**, 1853 (1994).

² W.H. Chan, R. Yuan, Ion-selective electrodes based on calix[4]arene tetraester in the determination of formaldehyde via in situ generation of ionic lipophilic hydrazone. **Analyst** **120**, 1055 (1995).