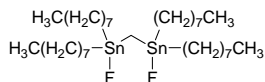


## Fluoride

### Fluoride Ionophore I

(*n*-Octyl)<sub>2</sub>(F)SnCH<sub>2</sub>Sn(F)(*n*-Octyl)<sub>2</sub>; Bis[fluoro(di-*n*-dioctyl)stannyl]methane; Stannane, methylenbis[fluoro(di-*n*-octyl)]

C<sub>33</sub>H<sub>70</sub>F<sub>2</sub>Sn<sub>2</sub> M<sub>r</sub> 742.33 [767355-17-9]



[02536](#) Selectophore<sup>®</sup>, function tested 100 mg

## Electrochemical Transduction

### Ion-Selective Electrodes

#### Application 1 and Sensor Type <sup>1,2</sup>

Assay of F<sup>-</sup> activity with solvent polymeric electrodes based on Fluoride Ionophore I. The major interfering ions are H<sub>2</sub>PO<sub>4</sub><sup>2-</sup> and SCN<sup>-</sup>.

#### Recommended Cell Assembly

Reference || sample solution || ion-selective membrane | 0.01 M KF ([47073](#)), buffered with 0.001M MES pH 5.5 ([69892](#)) | AgCl, Ag

#### Recommended Membrane Composition

1.5 wt% Fluoride Ionophore I ([02536](#))  
 65.3 wt% Bis(2-ethylhexyl) sebacate (DOS) ([84818](#))  
 33.2 wt% Poly(vinyl chloride) high molecular weight ([81392](#))

#### Electrode Characteristics and Function

Selectivity coefficients  $\log K_{F, X}^{\text{Pot}}$  as obtained by the separate solution method (0.01 M of the potassium salts).

$\log K_{F, \text{SCN}}^{\text{Pot}}$	0.1	$\log K_{F, \text{NO}_2}^{\text{Pot}}$	-2.0
$\log K_{F, \text{Br}}^{\text{Pot}}$	-2.5	$\log K_{F, \text{NO}_3}^{\text{Pot}}$	-3.3
$\log K_{F, \text{Cl}}^{\text{Pot}}$	-3.0	$\log K_{F, \text{H}_2\text{PO}_4}^{\text{Pot}}$	0.9

Slope of linear regression: 57-58 mV/decade

Nernstian electrode response ( $3 \times 10^{-5}$  to  $10^{-1}$  M KF)

Detection limit:  $\log a_{F^-} \sim -4.5$

<sup>1</sup> K. Perdikaki, I. Tsagkatakis, N. Chaniotakis, R. Altmann, K. Jurkschat, G. Reeske, Selective fluoride recognition and potentiometric properties of ion-selective electrodes based on bis(halodiphenylstannyl)alkanes, **Anal. Chim. Acta** 467, 197 (2002).

<sup>2</sup> N. Chaniotakis, K. Jurkschat, D. Müller, K. Perdikaki, G. Reeske, Bis[di-*n*-alkyl(fluoro)stannyl]methanes, (R<sub>2</sub>FSn)<sub>2</sub>CH<sub>2</sub> (R = *n*-octyl, *n*-dodecyl): Stable Fluoride-Selective Carriers, **Eur. J. Inorg. Chem.** 11, 2283 (2004).