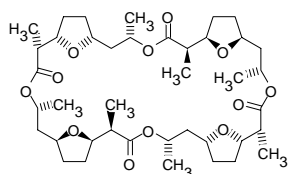


## Urea



### Ammonium ionophore I

(Nonactin\*)

$C_{40}H_{64}O_{12}$

$M_r$  736.95

[6833-84-7]

[09877](#)

**Selectophore<sup>®</sup>, function tested**

50 mg, 250 mg

\* purified and tested for the use in  $NH_4^+$ -selective electrodes

## Electrochemical Transduction

- Ion-Selective Electrodes

## Optical Transduction

## Electrochemical Transduction

### Ion-Selective Electrodes

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

Assay of  $\text{NH}_4^+$  activity derived from enzymatic decomposition of urea with solvent polymeric membrane electrodes based on Ammonium ionophore I.

Please note, this **Selectophore**<sup>®</sup> grade ionophore has been especially purified and tested for use in  $\text{NH}_4^+$ -selective electrodes.

#### Recommended Membrane Composition

0.60	wt%	Ammonium ionophore I ( <a href="#">09877</a> )
0.50	wt%	Tetraheptylammonium tetraphenylborate ( <a href="#">87293</a> )
63.60	wt%	Dibutyl sebacate ( <a href="#">84838</a> )
35.30	wt%	Poly(vinyl chloride) high molecular weight ( <a href="#">81392</a> )

#### Recommended Cell Assembly

Reference || sample solution || ion-selective membrane | 0.001 M  $\text{NH}_4\text{Cl}$  |  $\text{AgCl}$ ,  $\text{Ag}$

#### Electrode Characteristics and Function

Slope of linear regression: 53 mV ( $4 \cdot 10^{-6}$  to  $6 \cdot 10^{-3}$  M  $\text{NH}_4^+$  in 0.02 M TRIS/HCl,  $10^{-3}$  M  $\text{Na}_2\text{EDTA}$ ; pH 8)

Electrical resistance:  $\sim 3 \cdot 10^4 \Omega$

#### Application 2 and Sensor Type <sup>3</sup>

Assay of  $\text{NH}_4^+$  activity in an urea sensor which uses enzyme immobilisation on a nylon net together with a specially designed ammonium ion-selective field-effect transistor, coated with a solvent polymer membrane based on Ammonium ionophore I.

#### Recommended Membrane Composition

0.99	wt%	Ammonium ionophore I ( <a href="#">09877</a> )
0.60	wt%	Potassium tetrakis(4-chlorophenyl)borate ( <a href="#">60591</a> )
68.49	wt%	Tetraundecyl benzhydrol-3,3',4,4'-tetracarboxylate (ETH 2112) ( <a href="#">12103</a> )
29.82	wt%	Poly(vinyl chloride) high molecular weight ( <a href="#">81392</a> )

#### Recommended pH Buffer

0.001 M TRIS/HCl buffer; pH 7.3

#### Electrode Characteristics

Slope of linear regression: 57 mV ( $10^{-4}$  to  $3 \cdot 10^{-3}$  M urea)

Response time: 38s

Lifetime: >20d

#### Application 3 and Sensor Type <sup>4</sup>

Assay of  $\text{NH}_4^+$  activity derived from enzymatic decomposition of urea with solvent polymeric membrane electrodes based on Ammonium ionophore I.

#### Recommended Membrane Composition

4.00	wt%	Ammonium ionophore I ( <a href="#">09877</a> )
2.00	wt%	Palmitic acid ( <a href="#">76120</a> )
64.00	wt%	Bis(2-ethyl)hexyl sebacate (DOS) ( <a href="#">84818</a> )
30.00	wt%	Poly(vinyl chloride) high molecular weight ( <a href="#">81392</a> )

#### Recommended pH Buffer

Trizma base pH 7.00

#### Electrode Characteristics

Slope of linear regression:  $54.3 \pm 0.4$  mV ( $1.0 \cdot 10^{-5}$  to  $1.0 \cdot 10^{-2}$  M) urea;  $54.8 \pm 0.1$  mV ( $1.0 \cdot 10^{-6}$  to  $1.0 \cdot 10^{-1}$  M)  $\text{NH}_4^+$

Response time: < 1 min

Lifetime: >2 months

<sup>1</sup> S. Bilal Butt, K. Cammann, Enzyme Urea Biosensor Based on a Modified Potentiometric PVC-Nonactin Membrane Electrode for Assay of Urea in Blood. **Anal. Lett.** **25**, 1597 (1992).

<sup>2</sup> J.G. Schindler et al. Eur. Urea sensor for the continuous ammonium-selective enzymatic process control of the artificial kidney. **J. Clin. Chem. Clin. Biochem.** **32**, 145 (1994).

<sup>3</sup> S. Alegret, J. Bartolí, C. Jiménez, E. Martínez-Fabregas, D. Màrtorell, F. Valdés-Perezgasga, ISFET-based urea biosensor. **Sens. Actuators B15-16**, 453 (1993).

<sup>4</sup> E. Karakus, S. Pekyardimci, E. Kilic, Urea Biosensors Based on PVC Membrane Containing Palmitic Acid, **Artificial Cells, Blood Substitutes, and Biotechnology** **33**, 329 (2005).

## Optical Transduction

### Application 1 and Sensor Type<sup>5</sup>

Assay of  $\text{NH}_4^+$  activity in urea-sensitive reverse micelle based biooptode membranes. The ammonium-selective optode membrane contains Ammonium ionophore I and Chromoionophore III (ETH 5350).

#### Recommended Membrane Composition

0.37	wt%	Chromoionophore III ( <a href="#">27088</a> )
0.46	wt%	Ammonium ionophore I ( <a href="#">09877</a> )
0.62	wt%	Sodium tetrakis[3,5-bis(trifluoromethyl)phenyl]borate ( <a href="#">72017</a> )
5.76	wt%	Sodium bis(2-ethylhexyl)sulfosuccinate ( <a href="#">86139</a> )
0.49	wt%	Cyclohexanone ( <a href="#">29135</a> )
0.79	wt%	Urease ( <a href="#">94278</a> )
0.01	wt%	Water ( <a href="#">95283</a> )
58.70	wt%	Nitrophenyl octyl ether (NPOE) ( <a href="#">73732</a> )
32.80	wt%	Poly(vinyl chloride) high molecular weight ( <a href="#">81392</a> )

#### Recommended pH Buffer

0.018 M LiOH adjusted with phosphoric acid to pH 7.55

#### Optode Characteristics

Dynamic measuring range:  $10^{-5}$  to  $10^{-1}$  M urea at pH 7.55<sup>-</sup>

Response time: 95% response time 1-3.5 min

### Application 2 and Sensor Type<sup>6</sup>

Determination of urea based on the measurement of  $\text{NH}_4^+$  produced through enzymatic decomposition of urea, using fluorescence detection. The solvent polymeric optode membrane contains Ammonium ionophore I and Chromoionophore I.

#### Recommended Membrane Composition

9.50	wt%	Ammonium ionophore I ( <a href="#">09877</a> )
5.59	wt%	Chromoionophore I ( <a href="#">27086</a> )
4.47	wt%	Potassium tetrakis(4-chlorophenyl)borate ( <a href="#">60591</a> )
53.63	wt%	Bis(2-ethylhexyl)sebacate ( <a href="#">84818</a> )
26.81	wt%	Poly(vinyl chloride) high molecular weight ( <a href="#">81392</a> )

#### Recommended pH Buffer

10 or 200 mM TRIS/HCl buffer; pH 7.24.

#### Optode Characteristics

Dynamic range: 30  $\mu\text{M}$  to 10 mM  $\text{NH}_4^+$

Detection limit: 0.03 to 0.6 mM at near neutral pH

Response time: 4 min (enzyme membrane thickness 150  $\mu\text{m}$ )

### Application 3 and Sensor Type<sup>7</sup>

Determination of urease activity based on the measurement of  $\text{NH}_4^+$  produced through enzymatic decomposition of urea. This biosensor of urea incorporates an ammonium-selective bulk optode membrane containing Ammonium ionophore I and Chromoionophore III.

#### Recommended Membrane Composition

0.47	wt%	Ammonium ionophore I ( <a href="#">09877</a> )
0.43	wt%	Chromoionophore III ( <a href="#">27088</a> )
0.53	wt%	Sodium tetrakis[3,5-bis(trifluoromethyl)phenyl]borate ( <a href="#">72017</a> )
66.25	wt%	2-Nitrophenyl octyl ether ( <a href="#">73732</a> )
32.26	wt%	Poly(vinyl chloride) high molecular weight ( <a href="#">81392</a> )

#### Recommended pH Buffer

0.029 M NaOH/0.05 M  $\text{NaH}_2\text{PO}_4$ , pH 7.0.

#### Optode Characteristics

Dynamic range:  $10^{-4}$  to  $10^{-1}$  M  $\text{NH}_4\text{Cl}$  at pH 7.0

Detection limit: 0.03 to 0.6 mM at near neutral pH

Response time: 95% response time 0.5-2 min

<sup>5</sup> E. Vaillio, P. Walder, U.E. Spichiger, Development of micellar bioptodes membranes. **Anal. Meth. Instrum.** **2**, 145 (1995).

<sup>6</sup> O.S. Wolfbeis, H.Li, Fluorescence optical urea biosensor with an ammonium optrode as transducer. **Biosens. Bioelectron.** **8**, 161 (1993).

<sup>7</sup> C. Stamm, K. Seiler, W. Simon, Enzymatic biosensor for urea based on an ammonium ion-selective bulk optode membrane. **Anal. Chim. Acta** **282**, 239 (1993).