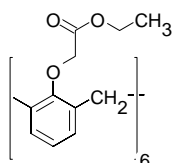


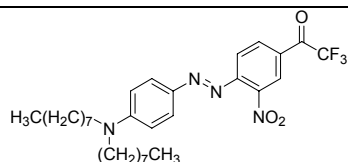
## Amine



### Amine ionophore I

(Calix[6]arene-hexaacetic acid hexaethyl ester)

 $C_{66}H_{72}O_{18}$      $M_r$  1153.28    [97600-45-8]

[06571](#)    **Selectophore®**, function tested    50 mg


### 4'-Diocetyl-amino-2-nitro-4-trifluoroacetylazobenzene

 (Chromoreactand CR-546; *N,N*-Diocetyl-amino-4'-trifluoroacetyl-2'-nitroazobenzene)

 $C_{30}H_{41}F_3N_4O_3$      $M_r$  562.67    [684281-90-1]

[08709](#)    **Selectophore®**    10 mg, 50 mg

## Electrochemical Transduction

### Ion-Selective Electrodes

### Optical Transduction

## Electrochemical Transduction

### Ion-Selective Electrodes

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

Determination of primary amine hydrochlorides by solvent polymeric membrane electrode.

#### Recommended Membrane Composition

5.00	wt%	Amine ionophore I ( <a href="#">06571</a> )
68.00	wt%	Bis(2-ethylhexyl) sebacate ( <a href="#">84818</a> )
27.00	wt%	Poly(vinyl chloride) high molecular weight ( <a href="#">81392</a> )

#### Recommended Cell Assembly

Reference || sample solution || ion-selective membrane | 0.01 M KCl | AgCl, Ag

#### Electrode Characteristics and Function

Selectivity coefficients  $\log K_{n - \text{Hexylammonium}, X}^{\text{Pot}}$  as obtained by the fixed interference method.

$\log K_{n - \text{Hexylammonium}, \text{Li}}^{\text{Pot}}$	-1.1
$\log K_{n - \text{Hexylammonium}, \text{Na}}^{\text{Pot}}$	-1.1
$\log K_{n - \text{Hexylammonium}, \text{K}}^{\text{Pot}}$	-1.1
$\log K_{n - \text{Hexylammonium}, \text{NH}_4}^{\text{Pot}}$	-1.0
$\log K_{n - \text{Hexylammonium}, \text{Butylammonium}}^{\text{Pot}}$	-1.0
$\log K_{n - \text{Hexylammonium}, \text{Piperidinium}}^{\text{Pot}}$	-1.1
$\log K_{n - \text{Hexylammonium}, \text{Diethylammonium}}^{\text{Pot}}$	-1.0
$\log K_{n - \text{Hexylammonium}, \text{Triethylammonium}}^{\text{Pot}}$	-1.4

Slope of linear regression: 57.0 mV ( $5 \cdot 10^{-5}$  –  $10^{-1}$  M n-hexylammonium chloride) (0.1 M TRIS-HCl buffer, pH 7).

Detection limit  $10^{-5}$  M n-hexylammonium chloride.

<sup>1</sup> W.H. Chan, K.K. Shiu, X.H. Gu, Ion-selective electrodes in organic analysis: primary amine selective polymeric membrane electrodes based on a calix[6]arene ionophore. **Analyst** **118**, 863 (1993).

<sup>2</sup> K. Odashima, K. Yagi, K. Tohda, Y. Umezawa, Potentiometric discrimination of organic amines by a liquid membrane electrode based on a lipophilic hexaester of calix[6]arene. **Anal. Chem.** **65**, 1074 (1993).

## Optical Transduction

### Application 1 and Sensor Type<sup>3</sup>

Determination of primary amines at pH 8 with solvent polymeric optode membranes based on Chromoionophore I (ETH 5294) and Amine ionophore I.

#### Recommended Membrane Composition

4.10	wt%	Amine ionophore I ( <a href="#">06571</a> )
0.47	wt%	Chromoionophore I ( <a href="#">27086</a> )
1.20	wt%	Sodium tetraphenylborate ( <a href="#">72018</a> )
62.80	wt%	Bis(2-ethylhexyl) phthalate ( <a href="#">80030</a> )
31.40	wt%	Poly(vinyl chloride) high molecular weight ( <a href="#">81392</a> )

#### Absorbance Maxima of Chromoionophore I in the membrane

$\lambda_{\text{max deprot.}}$	: 545 nm	$\lambda_{\text{max prot.}}$	: 600, 614 nm
--------------------------------	----------	------------------------------	---------------

#### Optode Characteristics and Function

Selectivity coefficients  $\log K_{n - \text{Octylammonium}, X}^{\text{Opt}}$  as obtained by the separate solution method (TRIS-HCl buffer pH 8).

$\log K_{n - \text{Octylammonium}, \text{Hexylammonium}}^{\text{Opt}}$	-1.15
$\log K_{n - \text{Octylammonium}, \text{Phenylethylammonium}}^{\text{Opt}}$	-1.30
$\log K_{n - \text{Octylammonium}, \text{Tetrabutylammonium}}^{\text{Opt}}$	-1.67
$\log K_{n - \text{Octylammonium}, n - \text{Butylammonium}}^{\text{Opt}}$	-1.95
$\log K_{n - \text{Octylammonium}, \text{Benzylammonium}}^{\text{Opt}}$	-2.10
$\log K_{n - \text{Octylammonium}, \text{Diethylammonium}}^{\text{Opt}}$	-3.53
$\log K_{n - \text{Octylammonium}, \text{Triethylammonium}}^{\text{Opt}}$	-3.30
$\log K_{n - \text{Octylammonium}, \text{NH}_4}^{\text{Opt}}$	-2.90
$\log K_{n - \text{Octylammonium}, \text{K}}^{\text{Opt}}$	-3.34
$\log K_{n - \text{Octylammonium}, \text{Na}}^{\text{Opt}}$	-4.90
$\log K_{n - \text{Octylammonium}, \text{Li}}^{\text{Opt}}$	-5.43

Detection range:  $10^{-6}$  –  $10^{-3}$  M octylamine at pH 8.0, down to  $10^{-7}$  at pH 8.5

Response time: 95% response time: 4 min ( $10^{-4}$  M to  $10^{-5}$  M)

Stability: After several hundred measurements the decrease in absorbance was 15%

<sup>3</sup> W.H. Chan, A.W.M. Lee, K. Wang, Design of a primary amine-selective optode membrane based on a lipophilic hexaester of calix[6]arene. **Analyst** **119**, 2809 (1994).

**Application 2 and Sensor Type**<sup>4</sup>

The chromoreactand 4'-Diocetyl-amino-2-nitro-4-trifluoroacetylazobenzene responds to dissolved amines provided a pH of above 10.0 is adjusted to obtain amines in the unprotonated form or when amines are present in the gaseous form.

**Recommended Membrane Composition**

0.83	wt%	4'-Diocetyl-amino-2-nitro-4-trifluoroacetylazobenzene ( <a href="#">08709</a> )
66.11	wt%	2-Nitrophenyl octyl ether ( <a href="#">73732</a> )
33.06	wt%	Poly(vinyl chloride) high molecular weight ( <a href="#">81392</a> )

**Preparation of the membrane**

1.0 mg of CR-546, 40 mg of PVC, and 80 mg of NPOE are dissolved in 0.7 ml of THF. 0.2 ml of this solution is pipetted on a rotating glass plate at 560 rpm to obtain layers of approximately 3-6  $\mu\text{m}$  thickness.

**Absorbance Maxima of 4'-Diocetyl-amino-2-nitro-4-trifluoroacetylazobenzene in the membrane**

$\lambda^{\text{max}}$  : 546 nm (in dichloromethane)

**Optode Characteristics and Function**

The response to, e.g. 1-propylamine, is in the 0.5 to 50 mM concentration range and the detection limit is 0.1 mM. Amines such as triethylamine, benzylamine or amphetamine can also be detected.

Response time: 2 min

<sup>4</sup> G. J. Mohr, Tailoring the sensitivity and spectral properties of a chromoreactand for the detection of amines and alcohols, **Anal. Chim. Acta** **508**, 233 (2004).