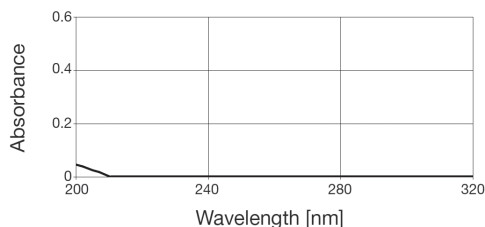


# LiChropur® ion-pair reagents

## What are ion-pair reagents?

These are ions with a strong hydrophobic character which form externally neutral associates with oppositely charged sample molecules. In this way it is possible to separate charged and uncharged molecules simultaneously.

The reagents in the LiChropur® range are specially selected for high UV transparency even at low detection wavelengths.

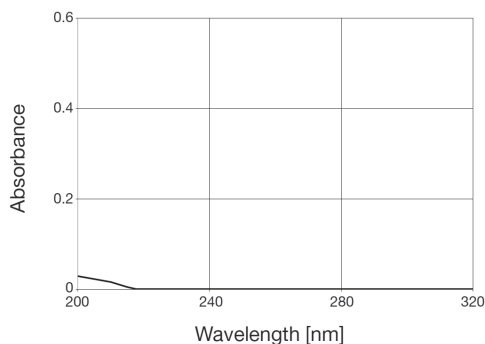


## UV spectrum of tetrabutylammonium hydrogen sulfate LiChropur®

Concentration: 0.005 mol/l in water

Reference: water

Path length: 1 cm



## UV spectrum of heptane-1-sulfonic acid sodium salt LiChropur®

Concentration: 0.005 mol/l in water

Reference: water

Path length: 1 cm

## With which columns and eluents can they be used?

They can be used fundamentally with all stationary phases: the eluent should contain at least 10% water as otherwise there is a risk (particularly with acetonitrile as the organic component) of precipitation.

When using long-chain ion-pair formers, e.g. cetyltrimethylammonium hydrogen sulfate or the sodium salt of dodecane sulfonic acid, the column to be used for the separation should be reserved purely for this purpose, as some irreversible adsorption to the stationary phase takes place, which causes changes in the separating performance.

## Instruction

For preparation of LiChropur® ion-pair reagents with buffer substance, we recommend the following instruction (according to the chromatographic requirement, the concentrations of LiChropur® ion-pair reagents can be optimized).

### Buffer substance pH 3.5:

#### 1. Composition:

50 mM  $\text{NaH}_2\text{PO}_4$ , 2.2 mM  $\text{NaHSO}_4$ ,  
5 mM ion-pair reagent [1-Butane sulfonic acid (Cat. No. 118303),  
pentane sulfonic acid (Cat. No. 1.18304),  
hexane sulfonic acid (Cat. No. 1.18305),  
heptane sulfonic acid (Cat. No. 1.18306), or  
octane sulfonic acid (Cat. No. 1.18307)]

or

50 mM  $\text{NaH}_2\text{PO}_4$ , 2.2 mM  $\text{NaHSO}_4$ ,  
0.5 mM ion-pair reagent [Dodecane sulfonic acid (Cat. No.1.18308)].

#### 2. Instruction:

Dissolve 138 g  $\text{NaH}_2\text{PO}_4 \times 1 \text{ H}_2\text{O}$  and  
6,07 g  $\text{NaHSO}_4 \times 1 \text{ H}_2\text{O}$  in  $\text{H}_2\text{O}$  and  
fill up to 1.0 l with  $\text{H}_2\text{O}$  = Solution 1.

Dissolve 5 mmol ion-pair reagent (1-Butane sulfonic acid, pentane sulfonic acid, hexane sulfonic acid, heptane sulfonic acid or octane sulfonic acid) in 50 ml of solution 1.  
Dilute with  $\text{H}_2\text{O}$  up to 1 l

or

dissolve 0.5 mmol ion-pair reagent (Dodecane sulfonic acid) in 50 ml of solution 1.  
Dilute with  $\text{H}_2\text{O}$  up to 1 l.

Buffer substance pH 6.5:

1. Composition:

50 mM Na<sub>2</sub>HPO<sub>4</sub>/NaH<sub>2</sub>PO<sub>4</sub> buffer +  
5 mM ion-pair reagent [Tetramethylammonium  
bisulfate (Cat. No. 87724, Sigma-Aldrich),  
Tetrabutylammonium hydrogen sulfate (Cat. No.  
1.18312)],

or

50 mM Na<sub>2</sub>HPO<sub>4</sub>/NaH<sub>2</sub>PO<sub>4</sub> buffer +  
0.5 mM ion-pair reagent [Cetyltrimethylammonium  
hydrogen sulfate (Cat. No. 52371, Sigma-  
Aldrich)].

Ordering information

Designation	Cat. No.	Content
1-Butanesulfonic acid sodium salt	1.18303.0025	25 g
	1.18303.0100	100 g
1-Pentanesulfonic acid sodium salt	1.18304.0025	25 g
	1.18304.0100	100 g
1-Hexanesulfonic acid sodium salt	1.18305.0025	25 g
	1.18305.0100	100 g
1-Heptanesulfonic acid sodium salt	1.18306.0025	25 g
	1.18306.0100	100 g
1-Octanesulfonic acid sodium salt	1.18307.0025	25 g
	1.18307.0100	100 g
	1.18307.1000	1000 g
1-Dodecanesulfonic acid sodium salt	1.18308.0025	25 g
	1.18308.0100	100 g
Tetramethylammonium bisulfate (Sigma-Aldrich)	87724-10G-F	10 g
	87724-50G-F	50 g
Tetrabutylammonium hydrogen sulfate	1.18312.0025	25 g
	1.18312.0100	100 g
Cetyltrimethylammonium hydrogen sulfate (Sigma-Aldrich)	52371-5G-F	5 g

Status: 2026-01-02

Made in Germany

Merck Life Science KGaA, 64271 Darmstadt, Germany,  
Tel. +49(0)6151 72-2440  
[www.sigmaaldrich.com/HPLC](http://www.sigmaaldrich.com/HPLC)



The life science business of Merck operates as MilliporeSigma in the U.S. and Canada.

© 2026 Merck KGaA, Darmstadt, Germany and/or its affiliates. All Rights Reserved. Merck, Supelco, Sigma-Aldrich, and LiChropur are trademarks of Merck KGaA, Darmstadt, Germany. All other trademarks are the property of their respective owners. Detailed information on trademarks is available via publicly available resources.