

Certificate

This certificate is designed in accordance with ISO Guide 31^[1]

Object of certification: **Sulfuric acid solution 0.1 mol/L in water**

Fluka Product No.: **68279 (Lot 42306A01)**

Composition: Sulfuric acid (puriss. p.a., 95-97%, Fluka Prod. No. 84720) dissolved in Ar-degassed high purity water (18.2 MΩ·cm, 0.2 μm filtered)

Certified value traceable to NIST and BAM certified reference materials and uncertainty according to ISO Guide 35 ^[2] and Eurachem/CITAC Guide ^[3]		
Constituent	Certified value at 20°C	Combined expanded uncertainty [$U = k u_c$; $k = 2$]
Sulfuric acid	100.0 mmol/L	0.4 mmol/L


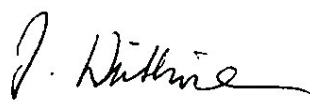

Intended use: Concentrate for preparation of eluents for ion chromatography

Storing and handling: This eluent concentrate solution shall be stored between 5°C and 30°C. In order to avoid evaporation the bottle should be tightly closed prior to use.

Expiry date: 28. May 2008 (unopened bottle)

Traceability statement: This eluent concentrate solution is traceable by potentiometric titration to NIST SRM 723c and also traceable to BAM certified titrimetric reference material (Fluka Prod. No. 93440).

Uncertainty calculation: All uncertainties are calculated according to Eurachem/CITAC Guide^[3] and reported as combined expanded uncertainties at the 95% confidence level. Contributions from reference material, potentiometric titration measurements and storing effects are included in the reported uncertainty budget.

Reviewing Chemist	Certification body	Date of release	Quality System
 K.-D. Schmidt, Ph.D.	 J. Wüthrich	June 15 th 2006	 SQS Reg. No. 16368-02

[1] ISO Guide 31, 1-7, 1st Ed. (1981), "Contents of certificates of reference materials"

[2] ISO Guide 35, 1-7, (2000), "Certification of reference materials – general and statistical principles"

[3] Eurachem/CITAC Guide, 1-120, 2nd Ed. (2000), "Quantifying uncertainty in analytical measurement"