



Specification

Certipur[®] Certified Reference Material

Manufacturer:	Merck KGaA, Frankfurter Str. 250, 64293 Darmstadt, Germany, Tel. +49(0)6151 720
Description of CRM:	Refractive index standard kit 3 1-Methylnaphthalene $n_D^{20} = 1.6160$ / Water $n_D^{20} = 1.3330$
Traceability:	This refractive index standard kit is measured by means of an Abbé-type refractometer and is directly traceable to the corresponding reference materials from NIST and PTB . <i>NIST: National Institute of Standards and Technology, Gaithersburg, USA</i> <i>PTB: Physikalisch Technische Bundesanstalt, Braunschweig, Germany</i>
Minimum shelf life:	3 years
Nominal value:	1-Methylnaphthalene $n_D^{20} = 1.6160$
Specification:	See table below

The certified value of the refractive index is measured lot specifically. Effective values can be found in the lot specific Certificate of Analysis. The measurement uncertainty is shown in the table below. Effective values of refractive index are always in the range shown as specification in the table below. Method of analysis is refractive index measured by an Abbé-type refractometer.

Art. No.	Substance	Specification n_D^{20}	Uncertainty n_D^{20}	Specification n_D^{25}	Uncertainty n_D^{25}
108963	1-Methylnaphthalene	1.61570 – 1.61750	± 0.00010	1.61345 – 1.61525	± 0.00010
	Water	1.33290 – 1.33308	± 0.00010	1.33234 – 1.33252	± 0.00010