

Marking

Characterization acc. ADR UN 1002, air, compressed, 2.2 Class 2, 1A

Essential properties

Colourless, odorless gas mixture, compressed, same weight as air

Cylinder Marking



shoulder:
light green

Symbols of Risks



gas, compressed

For additional safety information see Material-/safety data sheet No. *-N2-O2-0001A

Valves / Manifolds

Valve connection acc. to national standards

Recommended Manifolds Spectrolab FM 51, FM 52exact
Spectrocem FE 51 / FE 52exact



Specifications / Forms of delivery				
		5.0 HC-free	Scientific	
Composition				
N ₂	=	79,5	79,5	Vol. %
O ₂	=	20,5	20,5	Vol. %
Impurities				
H ₂ O	<	5	2	ppmv
THC (as CH ₄)	<	0,1	0,05	ppmv
CO ₂	<	0,5	0,1	ppmv
CO	<	-	0,1	ppmv
NO _x	<	0,1	0,01	ppmv
SO ₂	<	-	0,005	ppmv
H ₂ S	<	-	0,005	ppmv
Cylinders / Contents				
F 10 200 bar		2,0	-	m ³
F 50 200 bar		9,8	9,8	m ³
B 12* F 50 200 bar		117,1	-	m ³

Tolerance +/- 2,5 % rel.

Remarks

Gas chromatography: Zero gas, oxidizing gas for flame ionisation detectors (FID)