

Propene



also: Propylene

PURE GASES

Marking

CAS-Number 115-07-1

Characterization acc. ADR UN 1077, Propene, 2.1
Class 2, 2F

Cylinder Marking



shoulder:
red

Essential properties

Colourless, flammable gas, liquified, heavier than air

Symbols of Risks



highly flammable



gas, liquified

Physical Properties

molecular weight: 42,080 kg/kmol
gas density at 0°C and 1,013 bar: 1,9138 kg/m³
density ratio to air: 1,4802
vapour pressure at 20°C: 10,199 bar

For additional safety information see Material-/safety data sheet No. *-C3H6-105

Valves / Manifolds

Valve connection acc. to national standards

Recommended Manifolds Spectrolab control valve PN 40



Specifications / Cylinders

		2.5	
Composition			
C ₃ H ₆	>	99,5	Vol.-%
Impurities			
C ₃ H ₈	<	4000	ppmv
other HC	<	1000	ppmv
Cylinders / Contents			
F 27		11,0	kg
F 79		33,0	kg

Remarks

Applications:
Feedstock for polypropylene
Precursor for acetone, isopropanol, acrylonitrile and other organic compounds

Marking

CAS-Number 115-07-1

Characterization acc. ADR UN 1077, Propene, 2.1
Class 2, 2F

Cylinder Marking

shoulder:
red

Essential properties

Colourless, flammable gas, liquified, heavier than air

Symbols of Risks

highly flammable



gas, liquified

For additional safety information see Material-/safety data sheet No. *-C3H6-105

Description

Colourless, highly flammable, liquified gas. Forms explosive mixtures in air. Violent reactions with strong oxidizers. Catalytic polymerisation.

detection detector for flammable gases

Safety data

Explosion Range 1,8 - 11,2 Vol. %
Ignition Temperature 485 °C

Materials

Cylinders and valves: any usual materials
Seals: PTFE, PCTFE, PVDF, PP

Physical Properties	
molecular weight	42,080 kg/kmol
Critical Point	
temperature	365,57 K
Pressure	46,646 bar
density	0,22339 kg/l
Triple Point	
temperature	87,89 K
Pressure	9,5*10 ⁻⁹ bar
Boiling Point	
temperature	225,46 K; -47,7 °C
liquid density	0,60941 kg/l
evaporation heat	439,5 kJ/kg
vapour pressure at 20°C	10,199 bar
gas density at 0°C and 1,013 bar	1,9138 kg/m ³
density ratio to air	1,4802
gas density at 15°C and 1 bar	1,755 kg/m ³
Conversion Factor	
liquid at Ts to m ³ gas (15°C, 1 bar)	
Virial Coefficient	
Bn at 0°C	-18,7*10 ⁻³ bar ⁻¹
B30 at 30°C	-13,4*10 ⁻³ bar ⁻¹
Gaseous State at 25°C and 1 bar	
specific heat capacity cp	1,5306 kJ/kg K
thermal conductivity	170*10 ⁻⁴ W/m K
dynam. viscosity	8,5*10 ⁻⁶ Ns/m ²