

Chemical Safety Products

Choose from a wide range of Thermo Scientific Nalgene products specially designed for dispensing, labeling, disposing of hazardous waste, transporting and carrying chemical bottles. Cat. Nos. 2425 and 2436 comply with this U.S. OSHA Hazard Communication Standard, 29 CFR Part 1910.1200 (f) (4).*

The Anatomy of Nalgene® Right-To-Know Safety Products

Nalgene Right-To-Know safety products help your laboratory meet regulatory requirements for labeling chemical containers in the workplace. These symbols have been devised to provide instantly recognizable, standard symbols for the care, handling and storage of commonly-used laboratory chemicals. Nalgene Right-To-Know bottles and labeling incorporate the standard U.S. Department of Transportation (DOT), National Fire Protection Association (NFPA) and International Chemical Society (ICS) formulas. They also comply with U.S. Occupational Safety and Health Administration (OSHA) Hazardous Communication standards. Bottles are break-resistant, flexible, translucent and guaranteed leakproof. Labels are waterproof, chemical resistant and adhere well to plastic, glass and metal.

Chemical Name and Symbol

Chemical name in bold print (generic name in parentheses), followed by the ICS formula.

NFPA Diamond

Uses the standard NFPA codes, ranking hazards according to the chemical's reaction to the presence of fire. The red, blue and yellow diamonds use a rating scale of 4 to 0, with 4 representing the greatest hazard and 0 the least. The bottom diamond uses pictograms.

Top Diamond

- Red: Flash Point
- 4 - Below 70°F/21°C
- 3 - Below 100°F/38°C
- 2 - Below 200°F/93°C
- 1 - Above 200°F/93°C
- 0 - Non-flammable

Right-hand Diamond

- Yellow: Reactivity
- 4 - Explosive
- 3 - Shock and heat may detonate
- 2 - Violent change may occur
- 1 - Unstable if heated
- 0 - Normally stable

Left-hand Diamond

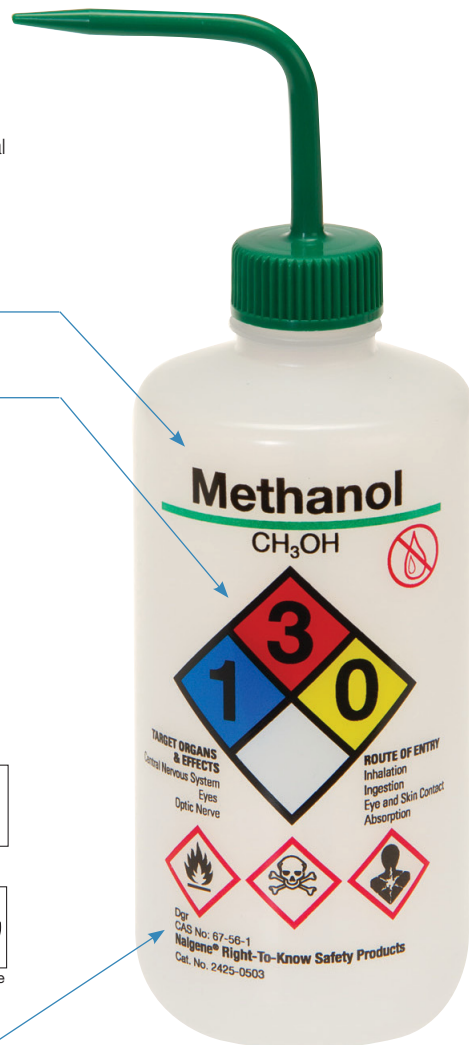
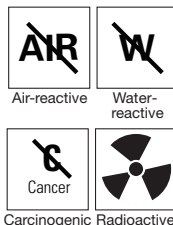
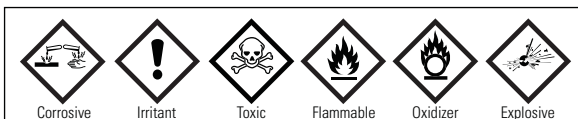
- Blue: Health Hazard
- 4 - Deadly
- 3 - Extremely dangerous
- 2 - Hazardous
- 1 - Slightly hazardous
- 0 - Normal material

Bottom Diamond

- White: Health Warnings
- Air-reactive
- Water-reactive
- Carcinogenic
- Radioactive

Hazard Code

Primary hazards represented by appropriate GHS symbol



DOT/UN Number

Four-digit number, preceded by UN or NA (North America), used to identify particular materials and regulate their transportation. These four digit numbers are called product identification numbers (PINs) under Canadian Transportation and Dangerous Goods Regulations.

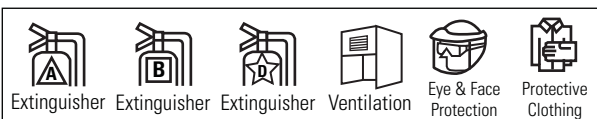
Hazard Code

(See description above)

Safety Equipment

Required protective equipment represented by the appropriate pictogram, including fire extinguishers:

- Extinguisher A - general use
- Extinguisher B - flammable liquids
- Extinguisher D - flammable materials



CAS Number

Chemical Abstract Service Number

Thermo Scientific Nalgene PolyPaper Right-to-Know Labels

Chemical: _____ DOT/UN No.: _____ CAS No.: _____

Hazard Codes: _____

Safety Equipment: _____

Prepared by: _____ Date: _____

Storage Color: Red Storage Color: White Storage Color: Yellow Storage Color: Orange Storage Color: Blue

Labels include: NFPA Code, Health Hazard, Flammability, Reactivity, Water-reactive, Air-reactive, Carcinogenic, Radioactive, and various GHS pictograms.

Storage Color

Color indicates the appropriate storage area:

- RED - flammable
- BLUE - health hazard
- YELLOW - reactive
- WHITE - contact hazard (corrosive)
- ORANGE - general storage

NFPA Diamond

(See description above)