

24-Hour Emergency Number
(VelocityEHS) 1-800-255-3924 in U.S.

cleaning@alconox.com

alconox.com

- [distributors](#)
- [technical information](#)
- [samples](#)
- [new developments](#)
- [white papers](#)
- [prices](#)

ALNOCHROMIX®

Oxidizing acid additive for glass cleaning

- Metal-free for trace metal glassware cleaning
- Surfactant-free for enzyme kinetics, cell and tissue culture cleaning
- Removes stubborn odors from containers
- Replaces carcinogenic and hazardous chromic acid
- Free rinsing to provide reliable results and no interfering residues

Used to clean: Laboratory glassware and other glass, and ceramic surfaces used for trace metal, enzyme kinetics, flavor profile, tissue and cell culture procedures..

Used to remove: Difficult organic, cyanide, and metal residues and odors

Surfaces cleaned: Glass, ceramic and acid resistant plastic surfaces.

Cleaning method: Soak or brush.

Directions: For making ALNOCHROMIX® solutions: Pour 1–2 pouches of ALNOCHROMIX® crystals into 1 large container (2.5L or 1gal capacity) of 98% concentrated sulfuric acid depending on desired cleaning strength. Stir lightly, crystals will dissolve in 10–15 minutes. If less sulfuric acid is used, use proportionally less ALNOCHROMIX®. Reseal tightly any unused crystals in a pouch. Store ALNOCHROMIX® sulfuric acid solution in a vented glass or HDPE containers.

For best results use freshly made solution within first 7 days of mixing. When cleaning solution is stale or discolored, it can be regenerated by adding new ALNOCHROMIX® cleaner.

Convenient Sizes	Alconox Cat. #
1 Box (10x3.1 Oz Packets)	2510
1 Case (4 Boxes)	2540
3 Cases (12 Boxes)	2512
Accessories:	
GI-45 Size Safety Vent Cap (3/Pk.)	2513

ALNOCHROMIX® cleaner is available from leading laboratory, hospital, clinical and industrial suppliers. To find a distributor for Alconox Inc. products, use the Locate Dealers search tool at alconox.com. To request samples, use the Order Samples form at alconox.com. Or call Alconox Inc. at +914.948.4040, today.

ALNTB.6

24-Hour Emergency Number
(VelocityEHS) 1-800-255-3924 in U.S.

cleaning@alconox.com

alconox.com

- distributors
- technical information
- samples
- new developments
- white papers
- prices

Physical Data

	Typical Value
Phosphate Content (as Phosphorus)	0%
Organic Carbon (% calculated w/w)	0%
Fragrance and Dye Content	0%
Color:	White
Form:	Powder
Shelf Life:	36 months from end of month of manufacture

Chemical Description

ALNOCHROMIX® cleaner consists primarily of inorganic persulfate oxidizing agent.

Important Warnings

- Do not close solution container airtight. **USE VENTED CAP** (vented caps cat no 2513).
- Never use nitric acid in place of sulfuric acid.
- Do not mix silver salt with ALNOCHROMIX®.
- Glassware containing silver salts should be rinsed before soaking in ALNOCHROMIX® and sulfuric acid solution.
- Wear suitable personal protective equipment for skin and eyes.

Health and Safety Information

RoHS Hazardous Ingredients:	None
EPA Priority Pollutants:	None
DOT Hazard Class:	UN1444 Ammonium Persulfate, 5.1, PGIII
Latex Content:	None in detergent, packaging materials or adhesives.
Eye Irritation:	Causes serious eye irritation. Wear appropriate protective equipment.
Skin Irritation:	Causes skin irritation. Wear appropriate protective equipment.
Carcinogenicity:	NTP = No IARC = No OSHA = No

All ingredients in ALNOCHROMIX® cleaner are listed in TSCA inventory.

Precautions

Contains inorganic persulfate. May cause allergy or asthma symptoms or breathing difficulties if inhaled. May cause respiratory irritation. Causes eye irritation. Causes skin irritation. May cause an allergic reaction. Harmful if swallowed. May intensify fire; oxidizer. Corrosive to metals and skin when mixed with sulfuric acid. A Safety Data Sheet is available at alconox.com. Contact Alconox Inc. for purchase specifications. Information presented is typical. Not to be taken as specifications. Typical data is not a specification.

While the information in this report should not be considered to be a product warranty, we urge you to investigate, test and verify the suitability of Alconox Inc. detergents for your specific application. We, of course, cannot give permission to use, or recommend the use of, our detergents where they infringe patents. No representation or warranty is made as to the safety of products or materials mentioned under the Federal Food Additives Amendment of 1958.

ALNTB.6