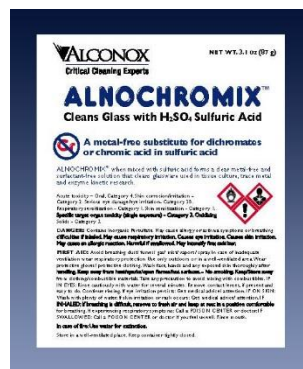


24-Hour Emergency Number
(ChemTel) 888.255.3924 in USA

cleaning@alconox.com

alconox.com

- Distributors
- Technical information
- Free samples
- New developments
- Whitepapers
- Prices



ALNOCHROMIX™

Cleans Glass with H₂SO₄ Sulfuric Acid

- 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 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™.

Catalog Number	Description
2510	1 BOX (10x3.1 oz PACKETS)
2540	1 CASE (4 BOXES)
2512	3 CASES (12 BOXES)
Accessories:	
2533	SAFETY VENT KIT (3 Silver safety gloves + 3 Vent Caps)
2513	SAFETY VENT CAP (3/pk.)
2521	1 PAIR SILVER SAFETY GLOVES

ALCONOX is available from leading laboratory, hospital, clinical and industrial suppliers. To find a distributor for Alconox, Inc. detergents, use the Locate Dealers search tool in the Support section of alconox.com. To request FREE samples, use the Get Samples form at alconox.com. Or call Alconox, Inc. at +914.948.4040, today. ALNTB.1

24-Hour Emergency Number
(ChemTel) 888.255.3924 in USA

cleaning@alconox.com

alconox.com

- Distributors
- Technical information
- Free samples
- New developments
- Whitepapers
- Prices

Physical Data

pH dilute in water (1%)	5
Specific gravity (g/ml)	1.0
Bulk Density (g/L.)	1000
Phosphate Content (as Phosphorus)	0%
Organic Carbon (% calculated w/w)	0%
Fragrance and Dye Content	0%
Color:	white
Form:	powder
Shelf Life:	2 years

Chemical Description

ALNOCHROMIX consists primarily of inorganic persulfate oxidizing agent

Important Warnings

- Do not close solution container airtight. Use vented cap (vented caps cat no 2513 and 2533).
- 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. Silver gloves Cat. No. 2521 and 2533 available.

Typical Data

Health 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 are listed in TSCA inventory.

Precautions

No special precautions other than good industrial hygiene and safety practices employed with any industrial chemical (see Directions). 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, can not 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.1