



Inhibitor-remover packings and ready-to-use, disposable prepacked columns offer a quick and convenient means of removing small amounts of inhibitors which are added to reagents or solvents that would otherwise be unstable (e.g., they may polymerize, oxidize or darken) on storage. The inhibitor removers are useful in applications which require that the stabilizer or inhibitor [i.e., hydroquinone (HQ), hydroquinone monomethyl ether (MEHQ, 4-methoxyphenol) or 4-*tert*-butylcatechol (TBC)] be removed prior to use.

TBC columns and packings are not recommended for use with polar solvents, including acrylic acid. We suggest vacuum distillation for such applications.

The prepacked glass columns come complete and ready for use. No prewash or pretreatment of any kind is necessary.

The TBC inhibitor removing column is moisture sensitive and comes packaged with end caps.

PROCEDURE

For optimum results, the following procedure is recommended.

STEP 1

Add the monomer (or solvent) to an addition funnel which is secured above the column.

STEP 2

Add the monomer **dropwise** to the column. Monitor the rate of addition to prevent overflow of the column. Collect the monomer in an appropriate container.

STEP 3

For low-melting solid monomers, keep the temperature of the column above the melting point during use. Do not overheat as this may cause the monomer to polymerize on the column. For very viscous or high-melting solid monomers, it is recommended that the monomer be diluted in an appropriate solvent before addition to the column.

STEP 4

Unused columns can be disposed of with other chemical solid waste. After use, the toxicity and hazards of the monomer from which the inhibitor is being removed, should be considered in deciding how to dispose of the packing material or column.

| INHIBITOR-REMOVER PRODUCTS | | | |
|-----------------------------------|--------------------------|--------------------|-------------------|
| Catalog No. | Inhibitor removed | Description | Dimensions |
| 30,631-2 | HQ, MEHQ | Column | 9 x 0.8 in. |
| 31,133-2 | HQ, MEHQ | Packing | |
| 30,632-0 | TBC | Column | 9 x 0.8 in. |
| 31,134-0 | TBC | Packing | |
| 34,019-7 | TBC | Column | 11 x 1.4 in. |



ALDRICH®

*Organics and Inorganics for
Chemical Synthesis*

1001 West Saint Paul Ave., Milwaukee, WI 53233

Telephone 414-273-3850 • 800-231-8327

Fax 414-273-4979 • 800-962-9591

Internet <http://www.sigma-aldrich.com>

E-mail aldrich@sial.com

© 1999 Sigma-Aldrich Co. Printed in the USA. Aldrich is a member of the Sigma-Aldrich Family. Sigma-Aldrich warrants that its products conform to the information contained in this and other Sigma-Aldrich publications. Purchaser must determine the suitability of the product for its particular use. See reverse side of invoice or packing slip for additional terms and conditions of sale.