

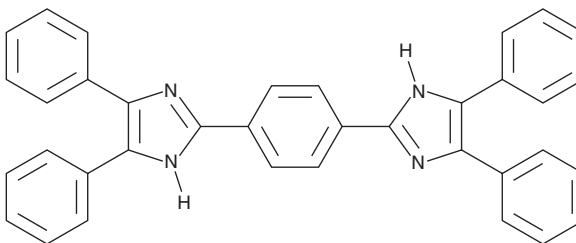
# PRODUCT INFORMATION



## MSX-130

Item No. 36804

CAS Registry No.: 4051-59-6  
Formal Name: 2,2'-(1,4-phenylene)bis[4,5-diphenyl-1H-imidazole]  
MF:  $C_{36}H_{26}N_4$   
FW: 514.6  
Purity:  $\geq 98\%$   
Supplied as: A solid  
Storage:  $-20^{\circ}\text{C}$   
Stability:  $\geq 4$  years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

MSX-130 is supplied as a solid. A stock solution may be made by dissolving the MSX-130 in the solvent of choice, which should be purged with an inert gas. MSX-130 is slightly soluble in DMSO and dimethyl formamide.

### Description

MSX-130 is an imidazole.<sup>1</sup> When used as the host in a host:guest complex with trichloroacetic acid, the complex displays excitation/emission maxima of 382/466 nm, respectively.

### Reference

1. Sakaino, Y., Shimizu, N., Takahashi, N., *et al.* Complex formation of imidazole derivatives and properties of the products. *Mol. Cryst. Liq. Cryst.* **187(1)**, 145-153 (1990).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 12/21/2022

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897  
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM