

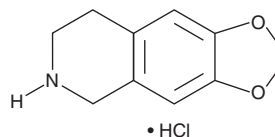
PRODUCT INFORMATION



TDIQ (hydrochloride)

Item No. 16083

CAS Registry No.: 15052-05-8
Formal Name: 5,6,7,8-tetrahydro-1,3-dioxolo[4,5-g]isoquinoline, monohydrochloride
Synonyms: MDTHIQ, 6,7-Methylenedioxy-1,2,3,4-tetrahydroisoquinoline
MF: $C_{10}H_{11}NO_2 \cdot HCl$
FW: 213.7
Purity: $\geq 95\%$
UV/Vis.: λ_{max} : 235, 292 nm
Supplied as: A crystalline solid
Stability: $-20^{\circ}C$
Storage: ≥ 5 years



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

Description

TDIQ (hydrochloride) is an analytical reference standard that is classified as a phenylalkylamine. It is related in structure to amphetamines but does not appear to affect locomotor activity.¹ TDIQ (hydrochloride) is a partial agonist of α_{2A} , α_{2B} , and α_{2C} -adrenergic receptors (K_i s = 75, 95, and 65 nM, respectively) but has low affinity for dopamine or serotonin receptors and does not affect catecholamine release in rodent brain synaptosomes.¹ Drug discrimination studies in rats indicate that it can serve as a discriminative stimulus, generalizing to cocaine (Item Nos. 16186, ISO60176) and partially to 3,4-MDMA (Item Nos. 13971, ISO60190) and methcathinone (Item No. 11709), but not to amphetamine (Item Nos 15797).¹ TDIQ has also been shown to exhibit anxiolytic and anorectic effects in rodent models.¹ This product is intended for forensic and research purposes.

Reference

1. Young, R. TDIQ (5,6,7,8-tetrahydro-1,3-dioxolo [4,5-g]isoquinoline): Discovery, pharmacological effects, and therapeutic potential. *CNS Drug Rev.* **13**(4), 405-422 (2007).

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, 03/03/2025

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