

Separation of stilbestrol derivatives

HPTLC Silica gel 60 RP-8 F254s

Plate	HPTLC pre-coated plate RP-8 254s F (1. 13725)
Eluent	Methanol/water 80/20 (v/v)
Migration Distance	5 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Diethylstilbestrol-dimethyl ether 2. Diethylstilbestrol-monomethyl ether 3. Diethylstilbestrol
Application volume	200 nl
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 254 nm

