

Separation of polystyrene 800

HPTLC Si 60 F254

Plate	TLC pre-coated plate gel 60 F254 with concentrating zone (P/N 1. 13728, alt P/N 1. 13727)
Eluent	Cyclohexane/toluene 75/25 (v/v)
Migration distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	Polystyrene standard 800
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 254 nm

