

Separation of insecticides

HPTLC Silica gel 60 F254

Plate	HPTLC pre-coated plate silica gel 60 F 254 (P/N: 1.05642, Alternative P/N: 1.05548)
Eluent	n-heptane
Migration Distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Endrin 2. DDT 3. Aldrin
Application volume	200 nl
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 254 nm

