

Separation of salicylic derivatives

TLC Silica gel 60 RP-18 F254s

Plate	TLC pre-coated plate RP-18 F 254s (P/N 1. 15389, alt. P/N 1. 15423)
Eluent	Methanol/IN acetic acid 50/50 (v/v)
Migration distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Salicylic acid 2. Acetylsalicylic acid 3. Salicyluric acid
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 254 nm

