

Separation of phenol carboxylic acids

HPTLC Si 60 F254

Plate	HPTLC pre-coated plate silica gel 60 F254 with concentrating zone (P/N 1. 13728, alt P/N 1. 13727)
Eluent	Acetonitrile/water 95/5 (v/v)
Migration distance	6 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Salicylic acid 0,15% 2. Syringic acid 0,05% 3. Ferulic acid 0,05% 4. o-Hydroxycinnamic acid 0,05%
Application volume	1 µl
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 265 nm

