

Separation of hydroxycinnamic 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/methanol/water 90/5/5 (v/v)
Migration distance	6 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. m-Hydroxycinnamic acid 2. p-Hydroxycinnamic acid 3. o-Hydroxycinnamic acid (all 0,02%)
Application volume	1 μ l
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 265 nm

