



1180 E. Ellsworth Rd, Ann Arbor, MI 48108
(800) 364-9897

Directions to Cayman Chemical:

Detroit Metro Airport

Take I-94 W
Take exit #177/State St
Continue on S State St
Turn left on Ellsworth Rd
Arrive at 1180 E Ellsworth Rd on the Right

Ann Arbor Municipal Airport

Turn Left onto State Rd
Turn Right onto Ellsworth
Arrive at 1180 E Ellsworth Rd on the Right

Detroit Willow Run Airport

Turn Left on Airport Dr
Turn Right on Tyler Rd
Take ramp toward Wiard Rd
Turn Right on Wiard Rd
Take Left ramp onto Ecorse Rd [US-12]
toward Ann Arbor
Take ramp onto I-94 West
Take exit #177/State St
Continue on S State St
Turn left on Ellsworth Rd
Arrive at 1180 E Ellsworth Rd on the Right

