


Coming From Delaware, Washington DC, Maryland and all points south:

1. Take I-95 North.
2. Exit 8 onto 202 North, Concord Pike (West Chester).
3. Go ~13 miles on 202 North and take a right side exit onto 202 North/322 West (Paoli and Downingtown).
4. Go 3 miles to the 322 West exit (Downingtown).
5. Go 3 miles (3 lights) to a stop sign and turn right continuing on 322 West.
6. Go 5 Miles to Boot Road (322 Truck West) and turn right.
7. Go approximately 0.5 miles on Boot Road. to the Ridgewood Place on the right.
8. Follow signs to Molecular Devices.

Coming From Philadelphia International Airport: ~35 miles

1. Take I-95 South (Delaware) from the airport.
2. Go 10 miles to exit 3A/322 West (West Chester).
3. Go 7 miles on 322 West to 1 South & 322 West (Baltimore Pike) and turn left.
4. Go one mile (four lights) and turn right on 322 West & 202 North.
5. Go 6 miles and take a right side exit onto 202 North/322 West (Paoli and Downingtown).
6. Go 3 miles to the 322 West exit (Downingtown).
7. Go 3 miles (3 lights) to a stop sign and turn right continuing on 322 West.
8. Go 5 Miles to Boot Road (322 Truck West) and turn right.
9. Go approximately 0.5 miles on Boot Road to Ridgewood Place on the right.
10. Follow signs to Molecular Devices.

Coming From Pennsylvania Turnpike:

1. Take PA Turnpike to exit 23/312 Downingtown.
2. Take Route 100 South for 1.5 miles, then right onto Route 113 South, Uwchlan Avenue.
3. Take 113 S for 4 miles. Continue past highway entrance to route 30, staying on 113 S.
4. When 113 S ends, take a right at the stop sign onto Business 30 West.
5. Almost immediately (200 feet) take a left onto Chestnut.
6. Stay on Chestnut, passing 2 stop signs and going up a hill and over a bridge.
7. When Chestnut ends at the traffic light, take a left onto Boot Road.
8. Take Boot Road for 1000 feet, past Skelp Level Road, then turn right into the Ridgewood Place.
9. Follow signs to Molecular Devices.