

From King of Prussia, PA:

Take Route 202 South.
Exit at Route 29 North (Great Valley).
Go through one light.
At 2nd light, make a Right onto Route 29 North.
At 2nd light, turn left onto Great Valley Parkway.
At the T junction, make a Left to stay on Great Valley Parkway.
We are on the Left at 271 Great Valley Parkway.

From New York or New Jersey:

Take the New Jersey Turnpike South or Route 1 South.
Exit onto the Pennsylvania Turnpike West.
Take the Pennsylvania Turnpike West to the Valley Forge Exit.
Take 202 South and follow directions from King of Prussia, PA.

From Washington DC and Baltimore:

Take 95 North.
As you approach Wilmington, DE keep to your left to stay on 95 North toward Philadelphia.
Soon after you reach Wilmington, Exit onto Route 202 North.
Follow Route 202 North for about 15 miles.
Exit onto Route 29 North (Great Valley).
At 2nd light, turn Left onto Great Valley Parkway.
At the T junction, make a Left to stay on Great Valley Parkway.
We are on the Left at 271 Great Valley Parkway.

From Philadelphia International Airport:

Follow route 95 South (Delaware), then route 476 North and then pick up route 76 West after 13 miles.
76 West will end, follow route 202 **South**, continue as instructed above from **King of Prussia**.

