

**EPIGENTEK SAME DAY DELIVERY RATE SHEET (as of January 2014)**

*SERVICE PROVIDED BY A1EXPRESS*

*Serving New York City and Long Island Areas*

**\$55**

11706, 11718, 11729, 11740, 11743, 11746, 11795

**\$65**

10474, 11001, 11003, 11004, 11005, 11010, 11020, 11021, 11022, 11023, 11024, 11030, 11040, 11042, 11050, 11096, 11101, 11102, 11103, 11104, 11105, 11106, 11222, 11351, 11352, 11354, 11355, 11356, 11357, 11358, 11359, 11360, 11361, 11362, 11363, 11364, 11365, 11366, 11367, 11368, 11369, 11370, 11371, 11372, 11373, 11374, 11375, 11377, 11378, 11379, 11380, 11381, 11385, 11386, 11405, 11411, 11412, 11413, 11415, 11416, 11417, 11418, 11419, 11420, 11421, 11422, 11423, 11425, 11426, 11427, 11428, 11429, 11430, 11431, 11432, 11433, 11434, 11435, 11436, 11439, 11451, 11501, 11507, 11509, 11510, 11514, 11516, 11518, 11520, 11530, 11531, 11535, 11542, 11545, 11547, 11548, 11549, 11550, 11551, 11552, 11553, 11554, 11555, 11556, 11557, 11558, 11559, 11560, 11561, 11563, 11565, 11566, 11568, 11569, 11570, 11571, 11572, 11575, 11576, 11577, 11579, 11580, 11581, 11590, 11592, 11594, 11595, 11596, 11597, 11598, 11599, 11691, 11692, 11693, 11694, 11697, 11701, 11703, 11704, 11707, 11708, 11709, 11710, 11714, 11717, 11721, 11724, 11726, 11731, 11732, 11735, 11736, 11737, 11747, 11750, 11751, 11753, 11756, 11757, 11758, 11762, 11765, 11771, 11773, 11774, 11775, 11783, 11791, 11793, 11797, 11798, 11801, 11802, 11803, 11804, 11815, 11854, 11855

**\$75**

11536, 11702, 11722, 11725, 11730, 11749, 11752, 11760, 11788

**\$85**

10001, 10002, 10003, 10009, 10010, 10011, 10012, 10013, 10016, 10017, 10018, 10019, 10020, 10021, 10022, 10023, 10024, 10025, 10026, 10027, 10028, 10029, 10030, 10031, 10032, 10033, 10034, 10035, 10036, 10037, 10038, 10039, 10040, 10055, 10065, 10069, 10075, 10080, 10082, 10095, 10101, 10102, 10103, 10104, 10105, 10106, 10107, 10108, 10109, 10110, 10111, 10112, 10117, 10118, 10119, 10120, 10121, 10122, 10123, 10128, 10129, 10138, 10149, 10156, 10157, 10158, 10159, 10160, 10165, 10166, 10168, 10169, 10170, 10172, 10173, 10174, 10175, 10176, 10178, 10179, 10185, 10276, 10451, 10452, 10453, 10454, 10455, 10456, 10457, 10458, 10459, 10460, 10461, 10462, 10463, 10464, 10465, 10466, 10467, 10468, 10469, 10470, 10471, 10472, 10473, 10475, 10502, 10503, 10522, 10523, 10528, 10530, 10532, 10533, 10538, 10543, 10550, 10552, 10553, 10557, 10558, 10573, 10577, 10580, 10583, 10591, 10595, 10601, 10602, 10603, 10604, 10605, 10606, 10607, 10610, 10701, 10703, 10704, 10705, 10706, 10707, 10708, 10709, 10710, 10801, 10803, 10804, 10805, 11041, 11201, 11203, 11204, 11205, 11206, 11207, 11208, 11210, 11211, 11212, 11213, 11215, 11216, 11217, 11218, 11221, 11223, 11225, 11226, 11229, 11230, 11233, 11234, 11235, 11236, 11237, 11238, 11239, 11241, 11242, 11243, 11245, 11247, 11251, 11414, 11582, 11716, 11739, 11754, 11769, 11779, 11782, 11787, 11796

**\$95**

00501, 00544, 11002, 11051, 11052, 11053, 11054, 11055, 11499, 11690, 11741, 11755, 11767, 11768, 11770, 11780

**\$105**

11705, 11715, 11720, 11738, 11742, 11763, 11772, 11790, 11794

**\$115**

10499, 11256, 11424, 11713, 11727, 11733, 11776, 11777, 11784

**\$125**

07010, 07020, 07024, 07605, 07620, 07626, 07631, 07632, 07670, 11695, 11719, 11764, 11766, 11789, 11953, 11967, 11980

**\$135**

10044, 10087, 10151, 10153, 10154, 10155, 10163, 10164, 10167, 10171, 10177, 10259, 10261, 10301, 10314, 10551, 11109, 11778, 11973

**\$145**

10043, 10072, 10113, 10114, 10115, 10116, 10124, 10125, 10126, 10130, 10131, 10132, 10133, 10150, 10152, 10162, 10197, 10199, 10802, 11202, 11249, 11786, 11951, 11961

**\$155**

10702

**\$185**

06836, 10964