

# Commonly Abused Drugs (Base)

Entry Num	Library ID	Name	Entry Num	Library ID	Name
288	AB-288	"PEACE PILL"			
4	AB-4	1,1-DIPHENYL-1-(2-DIMETHYL-LAMINOPROPYL)-2-BUTANONE HYDROCHLORIDE	20	AB-20	ENE)-17-METHYLANDROSTAN-3-ONE
			24	AB-24	2,2,2-TRICHLORO-1,1-ETHANEDIOL
			25	AB-25	2,2,2-TRICHLORO-1,1-ETHANEDIOL
6	AB-6	1,1-DIPHENYL-1-(2-DIMETHYL-LAMINOPROPYL)-2-BUTANONE HYDROCHLORIDE	261	AB-261	2,2,2-TRICHLORO-1,1-ETHANEDIOL
			109	AB-109	2,4-DIOXO-3,3-DIETHYL-5-METHYLPIPERIDINE
270	AB-270	1,2,3,4,5,6-HEXAHYDRO-6,11-DIMETHYL-3-(3-METHYL-2-BUTENYL)-2,6-METHANO-3-BENZAZOCIN-8-OL	110	AB-110	2,4-DIOXO-3,3-DIETHYL-5-METHYLPIPERIDINE
			111	AB-111	2,4-DIOXO-3,3-DIETHYL-5-METHYLPIPERIDINE
171	AB-171	1,3,7-TRIMETHYLXANTHINE	39	AB-39	2,5-DIMETHOXY-alpha-METHYLPHENETHYLAMINE
181	AB-181	1,3,7-TRIMETHYLXANTHINE			
203	AB-203	1-(3-ANILINOPROPYL)-4-PHENYLISONIPECOTIC ACID, ETHYL ESTER	39	AB-39	2,5-DIMETHOXYAMPHETAMINE
			96	AB-96	2-(4-CHLOROPHENYL)TETRAHYDRO-3-METHYL-4H-1,3-THIAZIN-4-ONE
208	AB-208	1-(3-CYANO-3,3-DIPHENYL-PROPYL)-4-PHENYLISONIPECOTIC ACID, ETHYL ESTER, HYDROCHLORIDE	115	AB-115	1,1-DIOXIDE
					2-(4-CHLOROPHENYL)TETRAHYDRO-3-METHYL-4H-1,3-THIAZIN-4-ONE
211	AB-211	1-(p-AMINOPHENETHYL)-4-PHENYLISONIPECOTIC ACID, ETHYL ESTER	51	AB-51	1,1-DIOXIDE
					2-[(2,3-DIMETHYLPHENYL)AMINO]BENZOIC ACID
81	AB-81	1-[(3,4-DIMETHOXYPHENYL)METHYL]-6,7-DIMETHOXYISOQUINOLINE HYDROCHLORIDE	224	AB-224	2-[2-[(4-CHLOROPHENYL)PHENYL-METHYL]-1-PIPERAZINYL]ETHOXY]ETHANOL, HYDROCHLORIDE
211	AB-211	1-[2-(4-AMINOPHENYL)ETHYL]-4-PHENYL-4-PIPERIDINECARBOXYLIC ACID, ETHYL ESTER	225	AB-225	2-[2-[4-[(4-CHLOROPHENYL)PHENYLMETHYL]-1-PIPERAZINYL]ETHOXY]ETHANOL, HYDROCHLORIDE
221	AB-221	1-[2-(4-AMINOPHENYL)ETHYL]-4-PHENYL-4-PIPERIDINECARBOXYLIC ACID, ETHYL ESTER, DIHYDROCHLORIDE	59	AB-59	2-CHLORO-N,N-DIMETHYL-10H-PHENOTHIAZINE-10-PROPANAMINE, HYDROCHLORIDE
249	AB-249	1-CARBOXY-N,N,N-TRIMETHYLMETHANAMINIUM HYDROXIDE INNER SALT COMPOUND WITH 2,2,2-TRICHLORO-1,1-ETHANEDIOL (1:1)	60	AB-60	2-CHLORO-N,N-DIMETHYL-10H-PHENOTHIAZINE-10-PROPANAMINE, HYDROCHLORIDE
26	AB-26	1-CHLORO-3-ETHYL-1-PENTEN-4-YL-3-OL	61	AB-61	2-CHLORO-N,N-DIMETHYL-10H-PHENOTHIAZINE-10-PROPANAMINE, HYDROCHLORIDE
199	AB-199	1-CHLORO-3-ETHYL-1-PENTEN-4-YL-3-OL	62	AB-62	2-CHLORO-N,N-DIMETHYL-10H-PHENOTHIAZINE-10-PROPANAMINE, HYDROCHLORIDE
275	AB-275	1-CHLORO-3-ETHYL-1-PENTEN-4-YL-3-OL	125	AB-125	2-ETHYL-2-PHENYLGLUTARIMIDE
202	AB-202	1-ETHYNYLCYCLOHEXANOL CARBAMATE	127	AB-127	2-METHYL-2-PROPYL-1,3-PROPANEDIOL DICARBAMATE
74	AB-74	10,11-DIHYDRO-N,N-DIMETHYL-5H-DIBENZ[b,f]AZEPINE-5-PROPANAMINE	130	AB-130	2-METHYL-2-PROPYL-1,3-PROPANEDIOL DICARBAMATE
			131	AB-131	2-METHYL-2-PROPYL-1,3-PROPANEDIOL DICARBAMATE
67	AB-67	10,11-DIHYDRO-N,N-DIMETHYL-5H-DIBENZ[b,f]AZEPINE-5-PROPANAMINE, HYDROCHLORIDE	133	AB-133	2-METHYL-2-PROPYL-1,3-PROPANEDIOL DICARBAMATE
10	AB-10	10,11-DIHYDRO-N,N-DIMETHYL-5H-DIBENZO[a,d]CYCLOHEPTENE-delta(SUP 5,gamma)PROPYLAMINE HYDROCHLORIDE	160	AB-160	2-METHYL-2-PROPYL-1,3-PROPANEDIOL DICARBAMATE
			86	AB-86	2-METHYL-3-(2-METHYLPHENYL)-4(3H)-QUINAZOLINONE
58	AB-58	10-[2-(1-METHYL-2-PIPERIDINYL)ETHYL]-2-(METHYLTHIO)-10H-PHENOTHIAZINE	87	AB-87	2-METHYL-3-(2-METHYLPHENYL)-4(3H)-QUINAZOLINONE
			109	AB-109	3,3-DIETHYL-5-METHYL-2,4-PIPERIDINEDIONE
227	AB-227	17-HYDROXY-2-(HYDROXYMETHYL-			

Entry Num	Library ID	Name	Entry Num	Library ID	Name
110	AB-110	3,3-DIETHYL-5-METHYL-2,4-PIPERIDINEDIONE	197	AB-197	5-ETHYL-1-METHYL-5-PHENYL-2,4,6(1H,3H,5H)-PYRIMIDINETRIONE
111	AB-111	3,3-DIETHYL-5-METHYL-2,4-PIPERIDINEDIONE	198	AB-198	5-ETHYL-1-METHYL-5-PHENYL-2,4,6(1H,3H,5H)-PYRIMIDINETRIONE
100	AB-100	3,7-DIHYDRO-1,3,7-TRIMETHYL-1H-PURINE-2,6-DIONE	128	AB-128	5-ETHYL-1-METHYL-5-PHENYL-2,4,6-(1H,3H,5H)-PYRIMIDINETRIONE
171	AB-171	3,7-DIHYDRO-1,3,7-TRIMETHYL-1H-PURINE-2,6-DIONE	129	AB-129	5-ETHYL-1-METHYL-5-PHENYL-2,4,6-(1H,3H,5H)-PYRIMIDINETRIONE
248	AB-248	3-(10,11-DIHYDRO-5H-DIBENZO[a,d]CYCLOHEPTEN-5-YLIDENE)-N-METHYL-1-PROPANAMINE HYDROCHLORIDE	168	AB-168	5-ETHYL-5-(1-METHYLBUTYL)-2,4,6(1H,3H,5H)-PYRIMIDINETRIONE MONOSODIUM SALT
125	AB-125	3-ETHYL-3-PHENYL-2,6-PIPERIDINEDIONE	169	AB-169	5-ETHYL-5-(1-METHYLBUTYL)-2,4,6(1H,3H,5H)-PYRIMIDINETRIONE MONOSODIUM SALT
157	AB-157	3-ETHYL-3-PHENYL-2,6-PIPERIDINEDIONE	134	AB-134	5-ETHYL-5-(1-METHYLBUTYL)-2,4,6(1H,3H,5H)PYRIMIDINETRIONE MONOSODIUM SALT
158	AB-158	3-ETHYL-3-PHENYL-2,6-PIPERIDINEDIONE	149	AB-149	5-ETHYL-5-(1-METHYLPROPYL)-2,4,6(1H,3H,5H)-PYRIMIDINETRIONE SODIUM SALT
159	AB-159	3-ETHYL-3-PHENYL-2,6-PIPERIDINEDIONE	150	AB-150	5-ETHYL-5-(1-METHYLPROPYL)-2,4,6(1H,3H,5H)-PYRIMIDINETRIONE SODIUM SALT
264	AB-264	3-METHYL-2-PHENYLMORPHOLINE HYDROCHLORIDE	151	AB-151	5-ETHYL-5-(1-METHYLPROPYL)-2,4,6(1H,3H,5H)-PYRIMIDINETRIONE SODIUM SALT
265	AB-265	3-METHYL-2-PHENYLMORPHOLINE HYDROCHLORIDE	153	AB-153	5-ETHYL-5-(1-METHYLPROPYL)-2,4,6(1H,3H,5H)-PYRIMIDINETRIONE SODIUM SALT
243	AB-243	4,5-EPOXY-3,14-DIHYDROXY-17-METHYLMORPHINAN-6-ONE HYDROCHLORIDE	141	AB-141	5-ETHYL-5-(3-METHYLBUTYL)-2,4,6-(1H,3H,5H)-PYRIMIDINETRIONE
230	AB-230	4,5-EPOXY-3-HYDROXY-17-METHYLMORPHINAN-6-ONE HYDROCHLORIDE	142	AB-142	5-ETHYL-5-(3-METHYLBUTYL)-2,4,6-(1H,3H,5H)-PYRIMIDINETRIONE
222	AB-222	4,5-EPOXY-3-HYDROXY-5,17-DIMETHYLMORPHINAN-6-ONE HYDROCHLORIDE	143	AB-143	5-ETHYL-5-(3-METHYLBUTYL)-2,4,6-(1H,3H,5H)-PYRIMIDINETRIONE
207	AB-207	4,5alpha-EPOXY-3-METHOXY-17-METHYLMORPHINAN-6-ONE TARTRATE (1:1)	144	AB-144	5-ETHYL-5-(3-METHYLBUTYL)-2,4,6-(1H,3H,5H)-PYRIMIDINETRIONE
83	AB-83	4,6-DIMETHYL-2-SULFANILAMINE DOPYRIMIDINE	186	AB-186	5-ETHYL-5-PHENYL-2,4,6(1H,3H,5H)-PYRIMIDINETRIONE MONOSODIUM SALT
8	AB-8	4-[2-(DIMETHYLAMINO)ETHYL]PHENOL	188	AB-188	5-ETHYL-5-PHENYL-2,4,6(1H,3H,5H)-PYRIMIDINETRIONE
164	AB-164	5-(1-CYCLOHEPTEN-1-YL)-5-ETHYL-2,4,6(1H,3H,5H)-PYRIMIDINETRIONE	189	AB-189	5-ETHYL-5-PHENYL-2,4,6(1H,3H,5H)-PYRIMIDINETRIONE
114	AB-114	5-(1-CYCLOHEXEN-1-YL)-1,5-DIMETHYL-2,4,6(1H,3H,5H)-PYRIMIDINETRIONE	196	AB-196	5-ETHYL-5-PHENYL-2,4,6(1H,3H,5H)-PYRIMIDINETRIONE
136	AB-136	5-(1-METHYLBUTYL)-5-(2-PROPENYL)-2,4,6(1H,3H,5H)-PYRIMIDINETRIONE MONOSODIUM SALT	257	AB-257	5-ETHYL-5-PHENYL-2,4,6(1H,3H,5H)-PYRIMIDINETRIONE
177	AB-177	5-(1-METHYLBUTYL)-5-(2-PROPENYL)-2,4,6(1H,3H,5H)-PYRIMIDINETRIONE MONOSODIUM SALT	279	AB-279	5-ETHYL-5-PHENYL-2,4,6(1H,3H,5H)-PYRIMIDINETRIONE
193	AB-193	5-(1-METHYLBUTYL)-5-(2-PROPENYL)-2,4,6(1H,3H,5H)-PYRIMIDINETRIONE MONOSODIUM SALT	188	AB-188	5-ETHYL-5-PHENYLBARBITURIC ACID
217	AB-217	5-(1-METHYLETHYL)-5-(2-PROPENYL)-2,4,6(1H,3H,5H)-PYRIMIDINETRIONE	189	AB-189	5-ETHYL-5-PHENYLBARBITURIC ACID
163	AB-163	5-(1-METHYLPROPYL)-5-(2-PROPENYL)-2,4,6(1H,3H,5H)-PYRIMIDINETRIONE	196	AB-196	5-ETHYL-5-PHENYLBARBITURIC ACID
26	AB-26	5-CHLORO-3-ETHYLPENT-1-YN-4-EN-3-OL	37	AB-37	7,8-DIDEHYDRO-4,5-EPOXY-17-
199	AB-199	5-CHLORO-3-ETHYLPENT-1-YN-4-EN-3-OL			
152	AB-152	5-ETHYKL-5-(1-METHYLPROPYL)-2,4,6(1H,3H,5H)-PYRIMIDINETRIONE SODIUM SALT			

Entry Num	Library ID	Name	Entry Num	Library ID	Name
43	AB-43	METHYLMORPHINAN-3,6-DIOL DIACETATE (ESTER)	283	AB-283	ARBITRARY MIXTURE: COCAINE BASE AND COCAINE HYDROCHLORIDE
17	AB-17	7,8-DIDEHYDRO-4,5-EPOXY-17-METHYLMORPHINAN-3,6-DIOL DIACETATE (ESTER) AND (8alpha,9R)-6'-METHOXY-CINCHONAN-9-OL, MONOHYDROCHLORIDE	282	AB-282	ARBITRARY MIXTURE: HEROIN AND PHENYLPROPANOLAMINE
120	AB-120	7,8-DIDEHYDRO-4,5-EPOXY-3-METHOXY-17-METHYL-(5alpha,6alpha)MORPHINAN-6-OL, PHOSPHATE (1:1) (SALT)	287	AB-287	ARBITRARY MIXTURE: HEROIN HYDROCHLORIDE AND CAFFEINE
140	AB-140	7-CHLORO-1,3-DIHYDRO-1-METHYL-5-PHENYL-2H-1,4-BENZODIAZEPIN-2-ONE	286	AB-286	ARBITRARY MIXTURE: HEROIN HYDROCHLORIDE, METHAMPHETAMINE HYDROCHLORIDE AND COCAINE HYDROCHLORIDE
276	AB-276	7-CHLORO-1,3-DIHYDRO-3-HYDROXY-5-PHENYL-2H-1,4-BENZODIAZEPIN-2-ONE	293	AB-293	ARBITRARY MIXTURE: HEROIN, QUININE HYDROCHLORIDE, METHAPYRILENE HYDROCHLORIDE AND MANNITOL
277	AB-277	7-CHLORO-1,3-DIHYDRO-3-HYDROXY-5-PHENYL-2H-1,4-BENZODIAZEPIN-2-ONE	297	AB-297	ARBITRARY MIXTURE: HEROIN, QUININE HYDROCHLORIDE, PROCAINE AND MANNITOL HYDROCHLORIDE
278	AB-278	7-CHLORO-1,3-DIHYDRO-3-HYDROXY-5-PHENYL-2H-1,4-BENZODIAZEPIN-2-ONE	290	AB-290	ARBITRARY MIXTURE: HEROIN, QUININE HYDROCHLORIDE, STRYCHNINE AND MANNITOL
89	AB-89	7-CHLORO-N-METHYL-5-PHENYL-3H-1,4-BENZODIAZEPIN-2-AMINE 4-OXIDE	298	AB-298	ARBITRARY MIXTURE: METHAMPHETAMINE AND ASCORBIC ACID
90	AB-90	7-CHLORO-N-METHYL-5-PHENYL-3H-1,4-BENZODIAZEPIN-2-AMINE 4-OXIDE	299	AB-299	ARBITRARY MIXTURE: METHAMPHETAMINE HYDROCHLORIDE AND AMPHETAMINE SULFATE
91	AB-91	7-CHLORO-N-METHYL-5-PHENYL-3H-1,4-BENZODIAZEPIN-2-AMINE 4-OXIDE	285	AB-285	ARBITRARY MIXTURE: METHAMPHETAMINE HYDROCHLORIDE AND COCAINE HYDROCHLORIDE
92	AB-92	7-CHLORO-N-METHYL-5-PHENYL-3H-1,4-BENZODIAZEPIN-2-AMINE 4-OXIDE	281	AB-281	ARBITRARY MIXTURE: METHAMPHETAMINE HYDROCHLORIDE AND METHENAMINE
88	AB-88	9,10-DIDEHYDRO-N,N-DIETHYL-6-METHYLERGOLINE-8beta-CARBOXAMIDE	284	AB-284	ARBITRARY MIXTURE: METHAMPHETAMINE HYDROCHLORIDE AND METHYLENEDIAMPHETAMINE HYDROCHLORIDE (MDA)
99	AB-99	A.P.C. WITH DEMROL	295	AB-295	ARBITRARY MIXTURE: METHAMPHETAMINE HYDROCHLORIDE AND PROCAINE HYDROCHLORIDE
37	AB-37	ACETOMORPHINE	296	AB-296	ARBITRARY MIXTURE: METHYLENEDIAMPHETAMINE HYDROCHLORIDE AND REDUCING SUGAR
100	AB-100	ALERT-PEP CAPSULES	300	AB-300	ARBITRARY MIXTURE: PHENCYCLIDINE AND PROCAINE HYDROCHLORIDE
231	AB-231	alpha-METHYLBENZENEETHANAMINE SULFATE	294	AB-294	ARBITRARY MIXTURE: SECOBARBITAL AND AMOBARBITAL
268	AB-268	alpha-PHENYL-2-PIPERIDINEACETIC ACID, METHYL ESTER, HYDROCHLORIDE	224	AB-224	ATARAX 10 MG.
217	AB-217	ALURATE 500 MG.	22	AB-22	ATARAX 100 MG.
203	AB-203	ALVODINE 50 MG.	225	AB-225	ATARAX 25 MG.
141	AB-141	AMOBARBITAL, U.S.P. 30 GM.	248	AB-248	AVENTYL HYDROCHLORIDE 10 MG.
144	AB-144	AMYTAL 1-1/2 GR.	9	AB-9	AVENTYL HYDROCHLORIDE 25 MG.
142	AB-142	AMYTAL 1/2 GR.	23	AB-23	B & O SUPRETTES 60/15
143	AB-143	AMYTAL 3/4 GR.	122	AB-122	BAMADEX
146	AB-146	AMYTAL SODIUM 0.2 GM.	249	AB-249	BETA-CHLOR
121	AB-121	AMYTAL SODIUM 1 GR.	28	AB-28	BIPHETAMINE 7-1/2
145	AB-145	AMYTAL SODIUM 65 MG.	123	AB-123	BIPHETAMINE-T-20
227	AB-227	ANADROL 2.5 MG.	228	AB-228	BONTRIL TIMED NO. 1
147	AB-147	APPETROL	229	AB-229	BONTRIL TIMED NO. 2
148	AB-148	APPETROL-S.R.	63	AB-63	BUFOTENINE BIOXALATE
289	AB-289	ARBITRARY MIXTURE: AMPHETAMINE SULFATE, BUTABARBITAL, PENTOBARBITAL, SECOBARBITAL AND PHENOBARBITAL	150	AB-150	BUTICAPS 3/4 GR.
292	AB-292	ARBITRARY MIXTURE: BUTABARBITAL AND SECOBARBITAL	149	AB-149	BUTISOL R-A 30 MG.
291	AB-291	ARBITRARY MIXTURE: BUTABARBITAL, PENTOBARBITAL AND SECOBARBITAL	153	AB-153	BUTISOL SODIUM 100 MG.
			151	AB-151	BUTISOL SODIUM 15 MG.
			152	AB-152	BUTISOL SODIUM 50 MG.
			181	AB-181	CAFFEINE
			250	AB-250	CALCIDRINE SYRUP
			260	AB-260	CANNABIS

Entry Num	Library ID	Name	Entry Num	Library ID	Name
155	AB-155	CARBRITAL ELIXIR	84	AB-84	EMPIRIN COMPOUND WITH CODEINE NO. 2
182	AB-182	CARBRITAL HALF STRENGTH P-D 376	104	AB-104	EMPIRIN COMPOUND WITH CODEINE NO. 3
154	AB-154	CARBRITAL P-D 372	93	AB-93	EMPIRIN COMPOUND WITH CODEINE NO. 4
251	AB-251	CHLOR-TRIMETON 12 MG.	183	AB-183	EPHEDRINE AND AMYTAL
53	AB-53	COCAINE & PROCAINE HYDROCHLORIDE	184	AB-184	EPHEDRINE AND AMYTAL
195	AB-195	COCAINE HYDROCHLORIDE	185	AB-185	EPHEDRINE AND SECONAL SODIUM
97	AB-97	CODALAN NO. 3	274	AB-274	EPHEDRINE SULFATE CAPSULES 25 MG.
17	AB-17	CODEINE PHOSPHATE 1 GR.	5	AB-5	EPHEDRINE SULFATE CAPSULES 50 MG.
69	AB-69	CODEINE SULFATE 30 MG.	126	AB-126	EQUAGESIC
18	AB-18	CODEINE SULFATE 60 MG.	127	AB-127	EQUANIL 400 MG.
12	AB-12	CODEMPIRAL NO. 2	160	AB-160	EQUANIL L-A
33	AB-33	CODEMPIRAL NO. 3	186	AB-186	ESKABARB SPANSULE 1 GR.
239	AB-239	COMPAZINE SPANSULES 30 MG.	257	AB-257	ESKABARB SPANSULE 1-1/2 GR.
252	AB-252	COMPAZINE SYRUP	161	AB-161	ESKAPHEN B
82	AB-82	COMPOZ	162	AB-162	ETHOBRAL
79	AB-79	COPAVIN	57	AB-57	ETHYLMORPHINE HYDROCHLORIDE USP
101	AB-101	CORICIDIN WITH CODEINE 1/2 GR.	94	AB-94	EXCEDRIN
49	AB-49	CORICIDIN WITH CODEINE 1/4 GR.	105	AB-105	FIORINAL
240	AB-240	COSANYL COUGH SYRUP	106	AB-106	FIORINAL
231	AB-231	D-alpha-METHYLPHENETHYLAMINE SULFATE	107	AB-107	FIORINAL WITH CODEINE NO. 2
1	AB-1	D-DESOXYEPHEDRINE HYDROCHLORIDE 1 GM.	187	AB-187	HARMONYL
246	AB-246	d-N,alpha-DIMETHYLPHENETHYLAMINE HYDROCHLORIDE	164	AB-164	HEPTADORM
117	AB-117	DALMANE 30 MG.	43	AB-43	HEROIN & QUININE HYDROCHLORIDE
13	AB-13	DAPRISAL	37	AB-37	HEROIN (FREE BASE)
34	AB-34	DARVO-TRAN	216	AB-216	HEXAHYDRO-1-METHYL-4-PHENYL-LAZEPINE-4-CARBOXYLIC ACID, ETHYL ESTER
218	AB-218	DARVON 32 MG.	8	AB-8	HORDENINE
204	AB-204	DARVON 65 MG.	209	AB-209	HYDROCORDONE BITARTRATE
14	AB-14	DARVON COMPOUND	258	AB-258	HYDRYLLIN COMPOUND
102	AB-102	DARVON COMPOUND 65	210	AB-210	HYPTRAN TABS
36	AB-36	DARVON WITH ASA	71	AB-71	IBOGAINE HYDROCHLORIDE
205	AB-205	DARVON-N 100 MG.	260	AB-260	INDIAN HEMP
108	AB-108	DASIN CAPSULES	259	AB-259	IONAMIN '30'
103	AB-103	DEMEROL APAP	31	AB-31	IONAMIN '15'
206	AB-206	DEMEROL HYDROCHLORIDE 100 MG.	24	AB-24	KESSODRATE CAPSULES 250 MG.
156	AB-156	DEPROL	20	AB-20	KESSODRATE CAPSULES 500 MG.
219	AB-219	DESOXYN 15 MG.	211	AB-211	LERITINE 25 MG.
220	AB-220	DESOXYN 5 MG.	221	AB-221	LERITINE DIHYDROCHLORIDE
124	AB-124	DEXAMYL	15	AB-15	LEVO-DROMORAN
253	AB-253	DEXAMYL SPANSULE NO. 1	269	AB-269	LEVOMOR
254	AB-254	DEXAMYL SPANSULE NO. 2	90	AB-90	LIBRITAB 10 MG.
255	AB-255	DEXEDRINE SPANSULE 15 MG.	89	AB-89	LIBRITAB 5 MG.
231	AB-231	DEXEDRINE SPANSULE 5 MG.	91	AB-91	LIBRITABS 25 MG.
37	AB-37	DIAMORPHINE	92	AB-92	LIBRIUM 25 MG.
2	AB-2	DIDREX 50 MG	3	AB-3	LOMOTIL
7	AB-7	DIETHYLTRYPTAMINE OXALATE	280	AB-280	LOMOTIL LIQUID
207	AB-207	DIHYDROCODEINON BITARTRATE, N.F.	163	AB-163	LOTUSATE 2 GR.
230	AB-230	DIHYDROMORPHINONE HYDROCHLORIDE, USP	88	AB-88	LSD
8	AB-8	DIMETHYLTRYPTAMINE	189	AB-189	LUMINAL 1-1/2 GR.
208	AB-208	DIPHENOXYLATE HYDROCHLORIDE	196	AB-196	LUMINAL 1/2 GR.
232	AB-232	DONNAGESIC NO. 1 EXTENTABS	188	AB-188	LUMINAL 1/4 GR.
241	AB-241	DONNAGESIC NO. 2 EXTENTABS	88	AB-88	LYSERGIDE
157	AB-157	DORIDEN 0.125 GM.	260	AB-260	MARIJUANA
158	AB-158	DORIDEN 0.25 GM.	129	AB-129	MEBARAL 1-1/2 GR.
125	AB-125	DORIDEN 0.5 GM.	197	AB-197	MEBARAL 1/2 GR.
159	AB-159	DORIDEN 0.5 GM.	198	AB-198	MEBARAL 3 GR.
40	AB-40	EDRISAL	128	AB-128	MEBARAL 50 MG.
50	AB-50	EDRISAL WITH CODEINE 1/2 GR.	164	AB-164	MEDOMIN 213 MG.
42	AB-42	EDRISAL WITH CODEINE 1/4 GR.	58	AB-58	MELLARIL 50 MG.
256	AB-256	ELAVIL HYDROCHLORIDE 10 MG.	213	AB-213	MEPERGAN FORTIS
10	AB-10	ELAVIL HYDROCHLORIDE 25 MG.	212	AB-212	MEPEGAN
11	AB-11	EMPIRAL	130	AB-130	MEPROSPAN-200 MG.

Entry Num	Library ID	Name	Entry Num	Library ID	Name
131	AB-131	MEPROSPAN-400 MG.	95	AB-95	PERCODAN
32	AB-32	MESCALINE ACID SULFATE	98	AB-98	PERCOGESIC-C
32	AB-32	MESCALINE SULFATE	16	AB-16	PHANTOS
6	AB-6	METHADONE HYDROCHLORIDE 10 MG.	45	AB-45	PHENAPHEN WITH CODEINE NO. 2
4	AB-4	METHADONE HYDROCHLORIDE 5 MG.	46	AB-46	PHENAPHEN WITH CODEINE NO. 3
242	AB-242	METHAMPHETAMINE HYDROCHLORIDE	47	AB-47	PHENAPHEN WITH CODEINE NO. 4
132	AB-132	METHAQUALONE HYDROCHLORIDE	263	AB-263	PHENAZOCINE HYDROBROMIDE 10 MG.
222	AB-222	METOPON 3 MG.	167	AB-167	PHENCYCLIDINE (PURE)
133	AB-133	MILTOWN 400 MG.	65	AB-65	PHENERGAN VC EXPECTORANT WITH CODEINE
233	AB-233	MORPHINE SULFATE PURE	279	AB-279	PHENOBARBITAL 1 GR.
246	AB-246	N,alpha-DIMETHYLBENZE-NEETHAMINE HYDROCHLORIDE	203	AB-203	PIMINODINE
219	AB-219	N,alpha-DIMETHYLBENZE-NEETHANAMINE HYDROCHLORIDE	199	AB-199	PLACIDYL 1 GM.
220	AB-220	N,alpha-DIMETHYLBENZE-NEETHANAMINE HYDROCHLORIDE	26	AB-26	PLACIDYL 100 MG.
242	AB-242	N,alpha-DIMETHYLBENZE-NEETHANAMINE HYDROCHLORIDE	275	AB-275	PLACIDYL 500 MG.
72	AB-72	N,N-DIMETHYL-N'-(PHENYL-METHYL)-N'-2-PYRIDINYL-1,2-ETHANEDIAMINE	76	AB-76	POLARAMINE SYRUP
234	AB-234	N,N-DIMETHYL-N'-(PHENYL-METHYL)-N'-2-PYRIDINYL-1,2-ETHANEDIAMINE	51	AB-51	PONSTEL 250 MG.
267	AB-267	N,N-DIMETHYL-N'-(PHENYL-METHYL)-N'-2-PYRIDINYL-1,2-ETHANEDIAMINE	245	AB-245	PRE-SATE 65 MG.
51	AB-51	N-(2,3-XYLYL)ANTHRANILIC ACID	264	AB-264	PRELUDIN ENDURETS 25 MG.
83	AB-83	N <sup>1</sup> -(4,6-DIMETHYL-2-PYRIDINYL)SULFANILAMIDE	265	AB-265	PRELUDIN ENDURETS 75 MG.
44	AB-44	NEBRALINE	175	AB-175	PROZINE CAPSULES
166	AB-166	NEMBU-DONNA 1/2	176	AB-176	PROZINE HALF STRENGTH CAPSULES
165	AB-165	NEMBU-DONNA 1/4	266	AB-266	PRYDON SPANSULE 0.4 MG.
190	AB-190	NEMBU-GESIC	214	AB-214	PRYDONNAL SPANSULE
85	AB-85	NEMBUDEINE 1/2 GR.	267	AB-267	PYRIBENZAMINE 100 MG.
134	AB-134	NEMBUTAL 1-1/2 GR.	72	AB-72	PYRIBENZAMINE 25 MG.
168	AB-168	NEMBUTAL 50 MG.	234	AB-234	PYRIBENZAMINE 50 MG.
169	AB-169	NEMBUTAL SODIUM CAPSULES 3/4 GR.	73	AB-73	PYRIBENZAMINE AND EPHEDRINE
191	AB-191	NEO-SEDAPHEN LIQUID	86	AB-86	QUAALUDE-150
170	AB-170	NIDAR	87	AB-87	QUAALUDE-300
171	AB-171	NO DOZ	215	AB-215	RITALIN 10 MG.
261	AB-261	NOCTEC 250 MG.	268	AB-268	RITALIN 20 MG.
25	AB-25	NOCTEC 500 MG.	21	AB-21	RITALIN 5 MG.
110	AB-110	NOLUDAR 200 MG.	77	AB-77	ROBITUSSIN A-C
111	AB-111	NOLUDAR 300 MG.	78	AB-78	ROBITUSSIN DM
109	AB-109	NOLUDAR 50 MG.	136	AB-136	SECOBARBITAL SODIUM 1-1/2 GR.
262	AB-262	NOVAHISTINE DH	177	AB-177	SECOBARBITAL SODIUM, U.S.P. 30 MG.
64	AB-64	NUCODAN	200	AB-200	SECONAL 3/4 GR.
243	AB-243	NUMORPHAN 10 MG.	193	AB-193	SECONAL SODIUM PULVULE 1-1/2 GR.
192	AB-192	OBEDRIN	277	AB-277	SERAX 10 MG.
172	AB-172	OBEDRIN-LA	278	AB-278	SERAX 15 MG.
173	AB-173	OBESTAT	276	AB-276	SERAX 30 MG.
54	AB-54	OPIUM, N.F.	80	AB-80	SINUTAB WITH CODEINE
75	AB-75	OXYCODONE HYDROCHLORIDE 50 MG.	66	AB-66	SLEEP-EZE
8	AB-8	p-HYDROXY-N,N-DIMETHYLPHENETHYLAMINE	193	AB-193	SODIUM 5-ALLYL-5-(1-METHYL-BUTYL)BARBITURATE
55	AB-55	PANTOPON 1/3 GR.	112	AB-112	SODIUM 5-ETHYL-5-(1-METHYL-BUTYL)BARBITURATE
81	AB-81	PAPAVERINE HYDROCHLORIDE 1-1/2 GR.	178	AB-178	SOMA COMPOUND
244	AB-244	PARACODIN	179	AB-179	SOMA COMPOUND WITH CODEINE
223	AB-223	PARAFON WITH CODEINE	114	AB-114	SOMBULEX
135	AB-135	PAREST-200	52	AB-52	SOMINEX
118	AB-118	PAREST-400	137	AB-137	SOMNAFAC CAPSULES
174	AB-174	PENTACINE	27	AB-27	STELAZINE 1 MG.
112	AB-112	PENTOBARBITAL SODIUM 1-1/2 GR.	56	AB-56	STELAZINE 10 MG.
113	AB-113	PERCOBARB	29	AB-29	STELAZINE 2 MG.
			30	AB-30	STELAZINE 5 MG.
			38	AB-38	STP IN METHANOL
			83	AB-83	SULFAMETHAZINE
			48	AB-48	SYNALGOS-DC CAPSULES
			19	AB-19	SYNATAN
			246	AB-246	SYNDROX TABLETS
			194	AB-194	SYNTIL
			238	AB-238	T.T CODEINE SULFATE, N.F. 15 MG. (1/4 GR.)
			70	AB-70	T.T CODEINE SULFATE, N.F. 30

Entry Num	Library ID	Name
		MG. (1/2 GR.)
270	AB-270	TALWIN 50 MG.
19	AB-19	TANPHETAMIN
119	AB-119	TENUATE DOSPAN 75 MG.
138	AB-138	TEPANIL 25 MG.
41	AB-41	THEBAINE ALKALOID 10 MG. (OPIUM DERIVATIVE)
59	AB-59	THORAZINE 10 MG.
62	AB-62	THORAZINE 100 MG.
60	AB-60	THORAZINE 25 MG.
61	AB-61	THORAZINE 50 MG.
67	AB-67	TOFRANIL 10 MG.
74	AB-74	TOFRANIL 25 MG.
68	AB-68	TOFRANIL 50 MG.
115	AB-115	TRANCOPAL 100 MG.
96	AB-96	TRANCOPAL 200 MG.
25	AB-25	TRICHLOROACETALDEHYDE MONOHY- DRATE
139	AB-139	TUINAL PULVULE 302
180	AB-180	TUINAL PULVULE 303
201	AB-201	TUINAL PULVULE 304
235	AB-235	TUSSEND
247	AB-247	TUSSIONEX
271	AB-271	TYLENOL WITH CODEINE NO. 2
272	AB-272	TYLENOL WITH CODEINE NO. 3
120	AB-120	VALIUM 10 MG.
140	AB-140	VALIUM 5 MG.
202	AB-202	VALMID 7-1/2 GR.
226	AB-226	VISTARIL 100 MG.
236	AB-236	VISTARIL 25 MG.
273	AB-273	VISTARIL 50 MG.
237	AB-237	WILPO 8 MG.
216	AB-216	ZACTANE 75 MG.
35	AB-35	ZACTIRIN AND ASPIRIN
116	AB-116	ZACTIRIN COMPOUND



**Bio-Rad  
Laboratories**

**Informatics Division**  
www.informatics.bio-rad.com

**U.S. Sales**

Phone: +1 888 5 BIO-RAD • E-mail: [informatics.usa@bio-rad.com](mailto:informatics.usa@bio-rad.com)

**Europe**

Phone: +44 20 8328 2555 • Free phone: 00800 78945000 • E-mail: [informatics.europe@bio-rad.com](mailto:informatics.europe@bio-rad.com)

**Japan**

Phone: +81 03 (5811) 6287 • E-mail: [informatics.nbr@j.bio-rad.com](mailto:informatics.nbr@j.bio-rad.com)

**Rest of World**

Phone: +1 215 382 7800 • E-mail: [informatics.row@bio-rad.com](mailto:informatics.row@bio-rad.com)