

# Interferon Gamma Monoclonal Mouse Antibody (G-30)

## Product Description

This antibody recognizes a protein of 20-25 kDa, identified as human interferon gamma (IFN- $\gamma$ ). Multiple forms exist due to variable glycosylation and under non-denaturing conditions due to dimers and tetramers. T lymphocytes and NK cells mainly produce IFN- $\gamma$ . It is a pleiotropic cytokine involved in the regulation of nearly all phases of immune and inflammatory responses, including the activation, growth and differentiation of T cell, B cells, macrophages, NK cells and other cell types such as endothelial cells and fibroblasts. It has weak anti-viral and anti-proliferative activity, and potentiates the antiviral and anti-tumor effects of IFN- $\gamma$  (type I interferon).

Primary antibodies are available purified, or with a selection of fluorescent CF® dyes and other labels. CF® dyes offer exceptional brightness and photostability. See the [CF® Dye Brochure](#) for more information. Note: Conjugates of blue fluorescent dyes like CF®405S and CF®405M are not recommended for detecting low abundance targets, because blue dyes have lower fluorescence and can give higher non-specific background than other dye colors.

**Stock status:** Because Biotium offers a large number of antibody and conjugation options, primary antibody conjugates may be made to order. Typical lead times are up to one week for CF® dye and biotin conjugates, and up to 2-3 weeks for fluorescent protein and enzyme conjugates. Please email [order@biotium.com](mailto:order@biotium.com) to inquire about stock status and lead times before placing your order.

### Product attributes

Antibody number	#0347
Antibody reactivity (target)	Interferon Gamma
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	G-30
Isotype	IgG1, kappa
Molecular weight	20-25 kDa
Synonyms	AMCF1; Beta thromboglobulin like protein; CXCL8; Chemokine ligand 8 (CXCL8); EMOX1; EMOX2; EMOX3; EMOX4; EMOX5; EMOX6; EMOX7; EMOX8; EMOX9; EMOX10; EMOX11; EMOX12; EMOX13; EMOX14; EMOX15; EMOX16; EMOX17; EMOX18; EMOX19; EMOX20; EMOX21; EMOX22; EMOX23; EMOX24; EMOX25; EMOX26; EMOX27; EMOX28; EMOX29; EMOX30; EMOX31; EMOX32; EMOX33; EMOX34; EMOX35; EMOX36; EMOX37; EMOX38; EMOX39; EMOX40; EMOX41; EMOX42; EMOX43; EMOX44; EMOX45; EMOX46; EMOX47; EMOX48; EMOX49; EMOX50; EMOX51; EMOX52; EMOX53; EMOX54; EMOX55; EMOX56; EMOX57; EMOX58; EMOX59; EMOX60; EMOX61; EMOX62; EMOX63; EMOX64; EMOX65; EMOX66; EMOX67; EMOX68; EMOX69; EMOX70; EMOX71; EMOX72; EMOX73; EMOX74; EMOX75; EMOX76; EMOX77; EMOX78; EMOX79; EMOX80; EMOX81; EMOX82; EMOX83; EMOX84; EMOX85; EMOX86; EMOX87; EMOX88; EMOX89; EMOX90; EMOX91; EMOX92; EMOX93; EMOX94; EMOX95; EMOX96; EMOX97; EMOX98; EMOX99; EMOX100; EMOX101; EMOX102; EMOX103; EMOX104; EMOX105; EMOX106; EMOX107; EMOX108; EMOX109; EMOX110; EMOX111; EMOX112; EMOX113; EMOX114; EMOX115; EMOX116; EMOX117; EMOX118; EMOX119; EMOX120; EMOX121; EMOX122; EMOX123; EMOX124; EMOX125; EMOX126; EMOX127; EMOX128; EMOX129; EMOX130; EMOX131; EMOX132; EMOX133; EMOX134; EMOX135; EMOX136; EMOX137; EMOX138; EMOX139; EMOX140; EMOX141; EMOX142; EMOX143; EMOX144; EMOX145; EMOX146; EMOX147; EMOX148; EMOX149; EMOX150; EMOX151; EMOX152; EMOX153; EMOX154; EMOX155; EMOX156; EMOX157; EMOX158; EMOX159; EMOX160; EMOX161; EMOX162; EMOX163; EMOX164; EMOX165; EMOX166; EMOX167; EMOX168; EMOX169; EMOX170; EMOX171; EMOX172; EMOX173; EMOX174; EMOX175; EMOX176; EMOX177; EMOX178; EMOX179; EMOX180; EMOX181; EMOX182; EMOX183; EMOX184; EMOX185; EMOX186; EMOX187; EMOX188; EMOX189; EMOX190; EMOX191; EMOX192; EMOX193; EMOX194; EMOX195; EMOX196; EMOX197; EMOX198; EMOX199; EMOX200; EMOX201; EMOX202; EMOX203; EMOX204; EMOX205; EMOX206; EMOX207; EMOX208; EMOX209; EMOX210; EMOX211; EMOX212; EMOX213; EMOX214; EMOX215; EMOX216; EMOX217; EMOX218; EMOX219; EMOX220; EMOX221; EMOX222; EMOX223; EMOX224; EMOX225; EMOX226; EMOX227; EMOX228; EMOX229; EMOX230; EMOX231; EMOX232; EMOX233; EMOX234; EMOX235; EMOX236; EMOX237; EMOX238; EMOX239; EMOX240; EMOX241; EMOX242; EMOX243; EMOX244; EMOX245; EMOX246; EMOX247; EMOX248; EMOX249; EMOX250; EMOX251; EMOX252; EMOX253; EMOX254; EMOX255; EMOX256; EMOX257; EMOX258; EMOX259; EMOX260; EMOX261; EMOX262; EMOX263; EMOX264; EMOX265; EMOX266; EMOX267; EMOX268; EMOX269; EMOX270; EMOX271; EMOX272; EMOX273; EMOX274; EMOX275; EMOX276; EMOX277; EMOX278; EMOX279; EMOX280; EMOX281; EMOX282; EMOX283; EMOX284; EMOX285; EMOX286; EMOX287; EMOX288; EMOX289; EMOX290; EMOX291; EMOX292; EMOX293; EMOX294; EMOX295; EMOX296; EMOX297; EMOX298; EMOX299; EMOX300; EMOX301; EMOX302; EMOX303; EMOX304; EMOX305; EMOX306; EMOX307; EMOX308; EMOX309; EMOX310; EMOX311; EMOX312; EMOX313; EMOX314; EMOX315; EMOX316; EMOX317; EMOX318; EMOX319; EMOX320; EMOX321; EMOX322; EMOX323; EMOX324; EMOX325; EMOX326; EMOX327; EMOX328; EMOX329; EMOX330; EMOX331; EMOX332; EMOX333; EMOX334; EMOX335; EMOX336; EMOX337; EMOX338; EMOX339; EMOX340; EMOX341; EMOX342; EMOX343; EMOX344; EMOX345; EMOX346; EMOX347; EMOX348; EMOX349; EMOX350; EMOX351; EMOX352; EMOX353; EMOX354; EMOX355; EMOX356; EMOX357; EMOX358; EMOX359; EMOX360; EMOX361; EMOX362; EMOX363; EMOX364; EMOX365; EMOX366; EMOX367; EMOX368; EMOX369; EMOX370; EMOX371; EMOX372; EMOX373; EMOX374; EMOX375; EMOX376; EMOX377; EMOX378; EMOX379; EMOX380; EMOX381; EMOX382; EMOX383; EMOX384; EMOX385; EMOX386; EMOX387; EMOX388; EMOX389; EMOX390; EMOX391; EMOX392; EMOX393; EMOX394; EMOX395; EMOX396; EMOX397; EMOX398; EMOX399; EMOX400; EMOX401; EMOX402; EMOX403; EMOX404; EMOX405; EMOX406; EMOX407; EMOX408; EMOX409; EMOX410; EMOX411; EMOX412; EMOX413; EMOX414; EMOX415; EMOX416; EMOX417; EMOX418; EMOX419; EMOX420; EMOX421; EMOX422; EMOX423; EMOX424; EMOX425; EMOX426; EMOX427; EMOX428; EMOX429; EMOX430; EMOX431; EMOX432; EMOX433; EMOX434; EMOX435; EMOX436; EMOX437; EMOX438; EMOX439; EMOX440; EMOX441; EMOX442; EMOX443; EMOX444; EMOX445; EMOX446; EMOX447; EMOX448; EMOX449; EMOX450; EMOX451; EMOX452; EMOX453; EMOX454; EMOX455; EMOX456; EMOX457; EMOX458; EMOX459; EMOX460; EMOX461; EMOX462; EMOX463; EMOX464; EMOX465; EMOX466; EMOX467; EMOX468; EMOX469; EMOX470; EMOX471; EMOX472; EMOX473; EMOX474; EMOX475; EMOX476; EMOX477; EMOX478; EMOX479; EMOX480; EMOX481; EMOX482; EMOX483; EMOX484; EMOX485; EMOX486; EMOX487; EMOX488; EMOX489; EMOX490; EMOX491; EMOX492; EMOX493; EMOX494; EMOX495; EMOX496; EMOX497; EMOX498; EMOX499; EMOX500; EMOX501; EMOX502; EMOX503; EMOX504; EMOX505; EMOX506; EMOX507; EMOX508; EMOX509; EMOX510; EMOX511; EMOX512; EMOX513; EMOX514; EMOX515; EMOX516; EMOX517; EMOX518; EMOX519; EMOX520; EMOX521; EMOX522; EMOX523; EMOX524; EMOX525; EMOX526; EMOX527; EMOX528; EMOX529; EMOX530; EMOX531; EMOX532; EMOX533; EMOX534; EMOX535; EMOX536; EMOX537; EMOX538; EMOX539; EMOX540; EMOX541; EMOX542; EMOX543; EMOX544; EMOX545; EMOX546; EMOX547; EMOX548; EMOX549; EMOX550; EMOX551; EMOX552; EMOX553; EMOX554; EMOX555; EMOX556; EMOX557; EMOX558; EMOX559; EMOX560; EMOX561; EMOX562; EMOX563; EMOX564; EMOX565; EMOX566; EMOX567; EMOX568; EMOX569; EMOX570; EMOX571; EMOX572; EMOX573; EMOX574; EMOX575; EMOX576; EMOX577; EMOX578; EMOX579; EMOX580; EMOX581; EMOX582; EMOX583; EMOX584; EMOX585; EMOX586; EMOX587; EMOX588; EMOX589; EMOX590; EMOX591; EMOX592; EMOX593; EMOX594; EMOX595; EMOX596; EMOX597; EMOX598; EMOX599; EMOX600; EMOX601; EMOX602; EMOX603; EMOX604; EMOX605; EMOX606; EMOX607; EMOX608; EMOX609; EMOX610; EMOX611; EMOX612; EMOX613; EMOX614; EMOX615; EMOX616; EMOX617; EMOX618; EMOX619; EMOX620; EMOX621; EMOX622; EMOX623; EMOX624; EMOX625; EMOX626; EMOX627; EMOX628; EMOX629; EMOX630; EMOX631; EMOX632; EMOX633; EMOX634; EMOX635; EMOX636; EMOX637; EMOX638; EMOX639; EMOX640; EMOX641; EMOX642; EMOX643; EMOX644; EMOX645; EMOX646; EMOX647; EMOX648; EMOX649; EMOX650; EMOX651; EMOX652; EMOX653; EMOX654; EMOX655; EMOX656; EMOX657; EMOX658; EMOX659; EMOX660; EMOX661; EMOX662; EMOX663; EMOX664; EMOX665; EMOX666; EMOX667; EMOX668; EMOX669; EMOX670; EMOX671; EMOX672; EMOX673; EMOX674; EMOX675; EMOX676; EMOX677; EMOX678; EMOX679; EMOX680; EMOX681; EMOX682; EMOX683; EMOX684; EMOX685; EMOX686; EMOX687; EMOX688; EMOX689; EMOX690; EMOX691; EMOX692; EMOX693; EMOX694; EMOX695; EMOX696; EMOX697; EMOX698; EMOX699; EMOX700; EMOX701; EMOX702; EMOX703; EMOX704; EMOX705; EMOX706; EMOX707; EMOX708; EMOX709; EMOX710; EMOX711; EMOX712; EMOX713; EMOX714; EMOX715; EMOX716; EMOX717; EMOX718; EMOX719; EMOX720; EMOX721; EMOX722; EMOX723; EMOX724; EMOX725; EMOX726; EMOX727; EMOX728; EMOX729; EMOX730; EMOX731; EMOX732; EMOX733; EMOX734; EMOX735; EMOX736; EMOX737; EMOX738; EMOX739; EMOX740; EMOX741; EMOX742; EMOX743; EMOX744; EMOX745; EMOX746; EMOX747; EMOX748; EMOX749; EMOX750; EMOX751; EMOX752; EMOX753; EMOX754; EMOX755; EMOX756; EMOX757; EMOX758; EMOX759; EMOX760; EMOX761; EMOX762; EMOX763; EMOX764; EMOX765; EMOX766; EMOX767; EMOX768; EMOX769; EMOX770; EMOX771; EMOX772; EMOX773; EMOX774; EMOX775; EMOX776; EMOX777; EMOX778; EMOX779; EMOX780; EMOX781; EMOX782; EMOX783; EMOX784; EMOX785; EMOX786; EMOX787; EMOX788; EMOX789; EMOX790; EMOX791; EMOX792; EMOX793; EMOX794; EMOX795; EMOX796; EMOX797; EMOX798; EMOX799; EMOX800; EMOX801; EMOX802; EMOX803; EMOX804; EMOX805; EMOX806; EMOX807; EMOX808; EMOX809; EMOX810; EMOX811; EMOX812; EMOX813; EMOX814; EMOX815; EMOX816; EMOX817; EMOX818; EMOX819; EMOX820; EMOX821; EMOX822; EMOX823; EMOX824; EMOX825; EMOX826; EMOX827; EMOX828; EMOX829; EMOX830; EMOX831; EMOX832; EMOX833; EMOX834; EMOX835; EMOX836; EMOX837; EMOX838; EMOX839; EMOX840; EMOX841; EMOX842; EMOX843; EMOX844; EMOX845; EMOX846; EMOX847; EMOX848; EMOX849; EMOX850; EMOX851; EMOX852; EMOX853; EMOX854; EMOX855; EMOX856; EMOX857; EMOX858; EMOX859; EMOX860; EMOX861; EMOX862; EMOX863; EMOX864; EMOX865; EMOX866; EMOX867; EMOX868; EMOX869; EMOX870; EMOX871; EMOX872; EMOX873; EMOX874; EMOX875; EMOX876; EMOX877; EMOX878; EMOX879; EMOX880; EMOX881; EMOX882; EMOX883; EMOX884; EMOX885; EMOX886; EMOX887; EMOX888; EMOX889; EMOX890; EMOX891; EMOX892; EMOX893; EMOX894; EMOX895; EMOX896; EMOX897; EMOX898; EMOX899; EMOX900; EMOX901; EMOX902; EMOX903; EMOX904; EMOX905; EMOX906; EMOX907; EMOX908; EMOX909; EMOX910; EMOX911; EMOX912; EMOX913; EMOX914; EMOX915; EMOX916; EMOX917; EMOX918; EMOX919; EMOX920; EMOX921; EMOX922; EMOX923; EMOX924; EMOX925; EMOX926; EMOX927; EMOX928; EMOX929; EMOX930; EMOX931; EMOX932; EMOX933; EMOX934; EMOX935; EMOX936; EMOX937; EMOX938; EMOX939; EMOX940; EMOX941; EMOX942; EMOX943; EMOX944; EMOX945; EMOX946; EMOX947; EMOX948; EMOX949; EMOX950; EMOX951; EMOX952; EMOX953; EMOX954; EMOX955; EMOX956; EMOX957; EMOX958; EMOX959; EMOX960; EMOX961; EMOX962; EMOX963; EMOX964; EMOX965; EMOX966; EMOX967; EMOX968; EMOX969; EMOX970; EMOX971; EMOX972; EMOX973; EMOX974; EMOX975; EMOX976; EMOX977; EMOX978; EMOX979; EMOX980; EMOX981; EMOX982; EMOX983; EMOX984; EMOX985; EMOX986; EMOX987; EMOX988; EMOX989; EMOX990; EMOX991; EMOX992; EMOX993; EMOX994; EMOX995; EMOX996; EMOX997; EMOX998; EMOX999; EMOX1000; EMOX1001; EMOX1002; EMOX1003; EMOX1004; EMOX1005; EMOX1006; EMOX1007; EMOX1008; EMOX1009; EMOX1010; EMOX1011; EMOX1012; EMOX1013; EMOX1014; EMOX1015; EMOX1016; EMOX1017; EMOX1018; EMOX1019; EMOX1020; EMOX1021; EMOX1022; EMOX1023; EMOX1024; EMOX1025; EMOX1026; EMOX1027; EMOX1028; EMOX1029; EMOX1030; EMOX1031; EMOX1032; EMOX1033; EMOX1034; EMOX1035; EMOX1036; EMOX1037; EMOX1038; EMOX1039; EMOX1040; EMOX1041; EMOX1042; EMOX1043; EMOX1044; EMOX1045; EMOX1046; EMOX1047; EMOX1048; EMOX1049; EMOX1050; EMOX1051; EMOX1052; EMOX1053; EMOX1054; EMOX1055; EMOX1056; EMOX1057; EMOX1058; EMOX1059; EMOX1060; EMOX1061; EMOX1062; EMOX1063; EMOX1064; EMOX1065; EMOX1066; EMOX1067; EMOX1068; EMOX1069; EMOX1070; EMOX1071; EMOX1072; EMOX1073; EMOX1074; EMOX1075; EMOX1076; EMOX1077; EMOX1078; EMOX1079; EMOX1080; EMOX1081; EMOX1082; EMOX1083; EMOX1084; EMOX1085; EMOX1086; EMOX1087; EMOX1088; EMOX1089; EMOX1090; EMOX1091; EMOX1092; EMOX1093; EMOX1094; EMOX1095; EMOX1096; EMOX1097; EMOX1098; EMOX1099; EMOX1100; EMOX1101; EMOX1102; EMOX1103; EMOX1104; EMOX1105; EMOX1106; EMOX1107; EMOX1108; EMOX1109; EMOX1110; EMOX1111; EMOX1112; EMOX1113; EMOX1114; EMOX1115; EMOX1116; EMOX1117; EMOX1118; EMOX1119; EMOX1120; EMOX1121; EMOX1122; EMOX1123; EMOX1124; EMOX1125; EMOX1126; EMOX1127; EMOX1128; EMOX1129; EMOX1130; EMOX1131; EMOX1132; EMOX1133; EMOX1134; EMOX1135; EMOX1136; EMOX1137; EMOX1138; EMOX1139; EMOX1140; EMOX1141; EMOX1142; EMOX1143; EMOX1144; EMOX1145; EMOX1146; EMOX1147; EMOX1148; EMOX1149; EMOX1150; EMOX1151; EMOX1152; EMOX1153; EMOX1154; EMOX1155; EMOX1156; EMOX1157; EMOX1158; EMOX1159; EMOX1160; EMOX1161; EMOX1162; EMOX1163; EMOX1164; EMOX1165; EMOX1166; EMOX1167; EMOX1168; EMOX1169; EMOX1170; EMOX1171; EMOX1172; EMOX1173; EMOX1174; EMOX1175; EMOX1176; EMOX1177; EMOX1178; EMOX1179; EMOX1180; EMOX1181; EMOX1182; EMOX1183; EMOX1184; EMOX1185; EMOX1186; EMOX1187; EMOX1188; EMOX1189; EMOX1190; EMOX1191; EMOX1192; EMOX1193; EMOX1194; EMOX1195; EMOX1196; EMOX1197; EMOX1198; EMOX1199; EMOX1200; EMOX1201; EMOX1202; EMOX1203; EMOX1204; EMOX1205; EMOX1206; EMOX1207; EMOX1208; EMOX1209; EMOX1210; EMOX1211; EMOX1212; EMOX1213; EMOX1214; EMOX1215; EMOX1216; EMOX1217; EMOX1218; EMOX1219; EMOX1220; EMOX1221; EMOX1222; EMOX1223; EMOX1224; EMOX1225; EMOX1226; EMOX1227; EMOX1228; EMOX1229; EMOX1230; EMOX1231; EMOX1232; EMOX1233; EMOX1234; EMOX1235; EMOX1236; EMOX1237; EMOX1238; EMOX1239; EMOX1240; EMOX1241; EMOX1242; EMOX1243; EMOX1244; EMOX1245; EMOX1246; EMOX1247; EMOX1248; EMOX1249; EMOX1250; EMOX1251; EMOX1252; EMOX1253; EMOX1254; EMOX1255; EMOX1256; EMOX1257; EMOX1258; EMOX1259; EMOX1260; EMOX1261; EMOX1262; EMOX1263; EMOX1264; EMOX1265; EMOX1266; EMOX1267; EMOX1268; EMOX1269; EMOX1270; EMOX1271; EMOX1272; EMOX1273; EMOX1274; EMOX1275; EMOX1276; EMOX1277; EMOX1278; EMOX1279; EMOX1280; EMOX1281; EMOX1282; EMOX1283; EMOX1284; EMOX1285; EMOX1286; EMOX1287; EMOX1288; EMOX1289; EMOX1290; EMOX1291; EMOX1292; EMOX1293; EMOX1294; EMOX1295; EMOX1296; EMOX1297; EMOX1298; EMOX1299; EMOX1300; EMOX1301; EMOX1302; EMOX1303; EMOX1304; EMOX1305; EMOX1306; EMOX1307; EMOX1308; EMOX1309; EMOX1310; EMOX1311; EMOX1312; EMOX1313; EMOX1314; EMOX1315; EMOX1316; EMOX1317; EMOX1318; EMOX1319; EMOX1320; EMOX1321; EMOX1322; EMOX1323; EMOX1324; EMOX1325; EMOX1326; EMOX1327; EMOX1328; EMOX1329; EMOX1330; EMOX1331; EMOX1332; EMOX1333; EMOX1334; EMOX1335; EMOX1336; EMOX1337; EMOX1338; EMOX1339; EMOX1340; EMOX1341; EMOX1342; EMOX1343; EMOX1344; EMOX1345; EMOX1346; EMOX1347; EMOX1348; EMOX1349; EMOX1350; EMOX1351; EMOX1352; EMOX1353; EMOX1354; EMOX1355; EMOX1356; EMOX1357; EMOX1358; EMOX1359; EMOX1360; EMOX1361; EMOX1362; EMOX1363; EMOX1364; EMOX1365; EMOX1366; EMOX1367; EMOX1368; EMOX1369; EMOX1370; EMOX1371; EMOX1372; EMOX1373; EMOX1374; EMOX1375; EMOX1376; EMOX1377; EMOX1378; EMOX1379; EMOX1380; EMOX1381; EMOX1382; EMOX1383; EMOX1384; EMOX1385; EMOX1386; EMOX1387; EMOX1388; EMOX1389; EMOX1390; EMOX1391; EMOX1392; EMOX1393; EMOX1394; EMOX1395; EMOX1396; EMOX1397; EMOX1398; EMOX1399; EMOX1400; EMOX1401; EMOX1402; EMOX1403; EMOX1404; EMOX1405; EMOX1406; EMOX1407; EMOX1408; EMOX1409; EMOX1410; EMOX1411; EMOX1412; EMOX1413; EMOX1414; EMOX1415; EMOX1416; EMOX1417; EMOX1418; EMOX1419; EMOX1420; EMOX1421; EMOX1422; EMOX1423; EMOX1424; EMOX1425; EMOX1426; EMOX1427; EMOX1428; EMOX1429; EMOX1430; EMOX1431; EMOX1432; EMOX1433; EMOX1434; EMOX1435; EMOX1436; EMOX1437; EMOX1438; EMOX1439; EMOX1440; EMOX1441; EMOX1442; EMOX1443; EMOX1444; EMOX1445; EMOX1446; EMOX1447; EMOX1448; EMOX1449; EMOX1450; EMOX1451; EMOX1452; EMOX1453; EMOX1454; EMOX1455; EMOX1456; EMOX1457; EMOX1458; EMOX1459; EMOX1460; EMOX1461; EMOX1462; EMOX1463; EMOX1464; EMOX1465; EMOX1466; EMOX1467; EMOX1468; EMOX1469; EMOX1470; EMOX1471; EMOX1472; EMOX1473; EMOX1474; EMOX1475; EMOX1476; EMOX1477; EMOX1478; EMOX1479; EMOX1480; EMOX1481; EMOX1482; EMOX1483; EMOX1484; EMOX1485; EMOX1486; EMOX1487; EMOX1488; EMOX1489; EMOX1490; EMOX1491; EMOX1492; EMOX1493; EMOX1494; EMOX1495; EMOX1496; EMOX1497; EMOX1498; EMOX1499; EMOX1500; EMOX1501; EMOX1502; EMOX1503; EMOX1504; EMOX1505; EMOX1506; EMOX1507; EMOX1508; EMOX1509; EMOX1510; EMOX1511; EMOX1512; EMOX1513; EMOX1514; EMOX1515; EMOX1516; EMOX1517; EMOX1518; EMOX1519; EMOX1520; EMOX1521; EMOX1522; EMOX1523; EMOX1524; EMOX1525; EMOX1526; EMOX1527; EMOX1528; EMOX1529; EMOX1530; EMOX1531; EMOX1532; EMOX1533; EMOX1534; EMOX1535; EMOX1536; EMOX1537; EMOX1538; EMOX1539; EMOX1540; EMOX1541; EMOX1542; EMOX1543; EMOX1544; EMOX1545; EMOX1546; EMOX1547; EMOX1548; EMOX1549;